

## Active Motif Epigenetic Services Publications

The papers below contain data generated by Active Motif's Epigenetic Services team. To learn more about our services, please give us a call or visit us at [www.activemotif.com/services](http://www.activemotif.com/services).

Technique	Target	Journal	Year	Reference
ChIP-Seq	RNA Pol II	Cell Reports	2022	Shin-Ichiro Ohno <i>et al.</i> Nuclear microRNAs release paused Pol II via the DDX21-CDK9 complex. <i>Cell Rep.</i> doi: 10.1016/j.celrep.2022.110673.
ChIP-Seq	H3K27ac	CRISPR J.	2022	Antonia A. Dominguez <i>et al.</i> CRISPR-Mediated Synergistic Epigenetic and Transcriptional Control. <i>CRISPR J.</i> doi: 10.1089/crispr.2021.0099.
ChIP-Seq	CTCF, H3K4me3, H3K27ac	Genome Research	2022	Blair W. Perry <i>et al.</i> Snake venom gene expression is coordinated by novel regulatory architecture and the integration of multiple co-opted vertebrate pathways. <i>Genome Res.</i> doi: 10.1101/gr.276251.121.
ChIP-Seq	SMARCA4, H3K-27me3	Cells	2022	Amelie Alfert <i>et al.</i> Smarcb1 Loss Results in a Deregulation of esBAF Binding and Impacts the Expression of Neurodevelopmental Genes. <i>Cells.</i> doi: 10.3390/cells11081354.
ChIP-Seq	BRD4	Biomed. Pharmacother.	2022	Laura M. Tsujikawa <i>et al.</i> Breaking boundaries: Pan BETi disrupt 3D chromatin structure, BD2-selective BETi are strictly epigenetic transcriptional regulators. <i>Biomed. Pharmacother.</i> doi: 10.1016/j.biopha.2022.113230.
ChIP-qPCR	H3K27ac, H3K-4me2	Nature Communications	2022	Emily V. Dutrow <i>et al.</i> Modeling uniquely human gene regulatory function via targeted humanization of the mouse genome. <i>Nat. Commun.</i> doi: 10.1038/s41467-021-27899-w.
ChIP-Seq	H3K4me3	Hepatol. Commun.	2022	Michael Schonfeld <i>et al.</i> Alcohol-associated fibrosis in females is mediated by female-specific activation of lysine demethylases KDM5B and KDM5C. <i>Hepatol. Commun.</i> doi: 10.1002/hep4.1967.
RIME	EOMES	iScience	2022	James A. Heslop <i>et al.</i> Chromatin remodeling is restricted by transient GATA6 binding during iPSC differentiation to definitive endoderm. <i>iScience.</i> doi: 10.1016/j.isci.2022.104300.
ChIP-Seq	H3K4me3	Hepatol. Commun.	2022	Michael Schonfeld <i>et al.</i> Male-Specific Activation of Lysine Demethylases 5B and 5C Mediates Alcohol-Induced Liver Injury and Hepatocyte Dedifferentiation. <i>Hepatol Commun.</i> doi: 10.1002/hep4.1895.
ChIP-Seq	BRD4, H3K27ac	Front Cell Dev Biol.	2022	Cho-Hao Lin <i>et al.</i> AZD5153, a Bivalent BRD4 Inhibitor, Suppresses Hepatocarcinogenesis by Altering BRD4 Chromosomal Landscape and Modulating the Transcriptome of HCC Cells. <i>Front Cell Dev Biol.</i> doi: 10.3389/fcell.2022.853652.
ATAC-Seq	—	Genes (Basel)	2022	Clarissa Boschiero <i>et al.</i> Characterization of Accessible Chromatin Regions in Cattle Rumen Epithelial Tissue during Weaning. <i>Genes (Basel)</i> doi: 10.3390/genes13030535.
ChIP-Seq	MAX	Cell Tissue Res.	2022	Weixu Ma <i>et al.</i> MAX deficiency impairs human endometrial decidualization through down-regulating OSR2 in women with recurrent spontaneous abortio. <i>Cell Tissue Res.</i> doi: 10.1007/s00441-022-03579-z.

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K4me1, H3K-4me3	FASEB J.	2022	Matthew Baxter <i>et al.</i> Circadian clock function does not require the histone methyltransferase MLL3. <i>FASEB J.</i> doi: 10.1096/fj.202200368R.
ATAC-Seq	—	Commun. Biol.	2022	Hidenori Machino <i>et al.</i> The metabolic stress-activated checkpoint LKB1-MARK3 axis acts as a tumor suppressor in high-grade serous ovarian carcinoma. <i>Commun. Biol.</i> doi: 10.1038/s42003-021-02992-4.
ChIP-Seq	H3K9me3	J. Immunol.	2022	Jun Hyung Sin <i>et al.</i> ATF7ip Targets Transposable Elements for H3K9me3 Deposition to Modify CD8+ T Cell Effector and Memory Responses. <i>J. Immunol.</i> doi: 10.4049/jimmunol.2100996.
ChIP-Seq	H3K9ac	Theranostics	2022	Grace Huang <i>et al.</i> Diabetes impairs cardioprotective function of endothelial progenitor cell-derived extracellular vesicles via H3K9Ac inhibition. <i>Theranostics.</i> doi: 10.7150/thno.70821.
ChIP-qPCR	WT1	Molecular Oncology	2022	Alain Chebly <i>et al.</i> Exploring hTERT promoter methylation in cutaneous T-cell lymphomas. <i>Mol Oncol.</i> doi: 10.1002/1878-0261.12946
ChIP-Seq	DNMT1	Gastroenterology	2022	Shikai Hu <i>et al.</i> NOTCH-YAP1/TEAD-DNMT1 Axis Drives Hepatocyte Reprogramming Into Intrahepatic Cholangiocarcinoma. <i>Gastroenterology.</i> doi: 10.1053/j.gastro.2022.05.007.
ATAC-Seq	—	Cell Rep. Med.	2022	Sally Mortlock <i>et al.</i> A multi-level investigation of the genetic relationship between endometriosis and ovarian cancer histotypes. <i>Cell Rep Med.</i> doi: 10.1016/j.xcrm.2022.100542.
ATAC-Seq, ChIP-Seq	H3K27ac, H3K-27me3, H3K4me1, H3K4me3, CTCF	Genomics	2022	Yahui Gae <i>et al.</i> Functional annotation of regulatory elements in cattle genome reveals the roles of extracellular interaction and dynamic change of chromatin states in rumen development during weaning. <i>Genomics.</i> doi: 10.1016/j.ygeno.2022.110296
ATAC-Seq, RNA-Seq	—	Autophagy	2022	Eutteum Jeong <i>et al.</i> The FACT complex facilitates expression of lysosomal and antioxidant genes through binding to TFEB and TFE3. <i>Autophagy.</i> doi: 10.1080/15548627.2022.2029671.
ChIP-Seq	H3K27ac	CRISPR J.	2022	Antonia A. Dominuez <i>et al.</i> CRISPR-mediated Synergistic Epigenetic and Transcriptional Control. <i>CRISPR J.</i> doi: 10.1089/crispr.2021.0099.
ChIP-Seq	YAP	J. Clin. Invest.	2022	Toshihide Kashihara <i>et al.</i> YAP mediates compensatory cardiac hypertrophy through aerobic glycolysis in response to pressure overload. <i>J. Clin. Invest.</i> doi: 10.1172/JCI150595.
ChIP-Seq	MAX	Cell Tissue Res.	2022	Weixu Ma <i>et al.</i> MAX deficiency impairs human endometrial decidualization through down-regulating OSR2 in women with recurrent spontaneous abortion. <i>Cell Tissue Res.</i> doi: 10.1007/s00441-022-03579-z.
ChIP-qPCR, Antibody Validation	KLF5	Science Advances	2022	Juliane Mooz <i>et al.</i> ARAF suppresses ERBB3 expression and metastasis in a subset of lung cancers. <i>Sci. Adv.</i> doi: 10.1126/sciadv.abk1538.
ChIP-Seq	H3K27ac, RNA pol II, ΔNp63	Nature Communications	2022	Marco Napoli <i>et al.</i> ΔNp63 regulates a common landscape of enhancer associated genes in non-small cell lung cancer. <i>Nat. Commun.</i> doi: 10.1038/s41467-022-28202-1.

Technique	Target	Journal	Year	Reference
ChIP-qPCR, Antibody Validation	H3K27ac, H3K-4me2	Nature Communications	2022	Emily V. Dutrow <i>et al.</i> Modeling uniquely human gene regulatory function via targeted humanization of the mouse genome. <i>Nat. Commun.</i> doi: 10.1038/s41467-021-27899-w.
ChIP-Seq	H3K4me3	Hepatol. Commun.	2022	Michael Schonfeld <i>et al.</i> Male-Specific Activation of Lysine Demethylases 5B and 5C Mediates Alcohol-Induced Liver Injury and Hepatocyte Dedifferentiation. <i>Hepatol. Commun.</i> doi: 10.1002/hep4.1895.
ATAC-Seq	—	Genes	2022	Clarissa Boschiero <i>et al.</i> Characterization of Accessible Chromatin Regions in Cattle Rumen Epithelial Tissue during Weaning. <i>Genes (Basel)</i> . doi: 10.3390/genes13030535.
ATAC-Seq	—	Mol. Cancer Res.	2022	Aditya Wirawan <i>et al.</i> A Novel Therapeutic Strategy Targeting the Mesenchymal Phenotype of Malignant Pleural Mesothelioma by Suppressing LSD1. <i>Mol. Cancer Res.</i> doi: 10.1158/1541-7786.MCR-21-0230.
ChIP-Seq	H3K9me3	J. Immunol.	2022	Jun Hyung Sin <i>et al.</i> ATF7ip Targets Transposable Elements for H3K9me3 Deposition to Modify CD8 + T Cell Effector and Memory Responses. <i>J. Immunol.</i> doi: 10.4049/jimmunol.2100996.
ATAC-Seq	—	Commun. Biol.	2022	Hidenori Machino <i>et al.</i> The metabolic stress-activated checkpoint LKB1-MARK3 axis acts as a tumor suppressor in high-grade serous ovarian carcinoma. <i>Commun. Biol.</i> doi: 10.1038/s42003-021-02992-4.
ATAC-Seq	—	Cancer Immunol. Res.	2022	Andrew Fedoriw <i>et al.</i> Inhibiting Type I arginine methyltransferase activity promotes the T cell mediated antitumor immune response. <i>Cancer Immunol. Res.</i> doi: 10.1158/2326-6066.CIR-21-0614.
MeDIP-Seq	—	EMBO Reports	2022	Antonia Piazzesi <i>et al.</i> CEST-2.2 overexpression alters lipid metabolism and extends longevity of mitochondrial mutants. <i>EMBO Rep.</i> doi: 10.15252/embr.202152606.
ChIP-qPCR	H3K27ac	RNA	2021	Ryan L. Setten <i>et al.</i> CRED9: A differentially expressed elncRNA regulates expression of transcription factor CEBPA. <i>RNA</i> . doi: 10.1261/rna.078752.121.
ATAC-Seq	—	EMBO Reports	2021	Raquel Sales-Gil <i>et al.</i> Non-redundant functions of H2A.Z.1 and H2A.Z.2 in chromosome segregation and cell cycle progression. <i>EMBO Rep.</i> doi: 10.15252/embr.202052061.
ATAC-Seq	—	Cell	2021	Florian Wimmers <i>et al.</i> The single-cell epigenomic and transcriptional landscape of immunity to influenza vaccination. <i>Cell</i> . doi: 10.1016/j.cell.2021.05.039.
Mod Spec	—	Blood	2021	Zachary C. Murphy <i>et al.</i> Regulation of RNA polymerase II activity is essential for terminal erythroid maturation. <i>Blood</i> . doi: 10.1182/blood.2020009903.
ATAC-Seq	—	J. Mol. Cell Cardiol.	2021	Douglas J. Chapski <i>et al.</i> Early adaptive chromatin remodeling events precede pathologic phenotypes and are reinforced in the failing heart. <i>J. Mol. Cell Cardiol.</i> doi: 10.1016/j.jmcc.2021.07.002.
ATAC-Seq	—	Nature Communications	2021	Joana Esteves de Lima <i>et al.</i> HIRA stabilizes skeletal muscle lineage identity. <i>Nat. Commun.</i> doi: 10.1038/s41467-021-23775-9.
ChIP-qPCR	LXRalpha	Br. J. Pharmacol.	2021	Louise Menegaut <i>et al.</i> Regulation of glycolytic genes in human macrophages by oxysterols: a potential role for liver X receptors. <i>Br. J. Pharmacol.</i> doi: 10.1111/bph.15358.

Technique	Target	Journal	Year	Reference
ChIP-Seq	Sp6	Biochem. Biophys. Res. Commun.	2021	Craig S. Rhodes <i>et al.</i> Sp6/Epiprofin is a master regulator in the developing tooth. <i>Biochem. Biophys. Res. Commun.</i> doi: 10.1016/j.bbrc.2021.10.017.
ChIP-Seq	BMAL1	Nature Communications	2021	Eleonore Maury <i>et al.</i> Circadian clock dysfunction in human omental fat links obesity to metabolic inflammation. <i>Nat. Commun.</i> doi: 10.1038/s41467-021-22571-9.
ChIP-Seq	p300	Clin. Transl. Med.	2021	Chuanjie Zhang <i>et al.</i> Epigenome screening highlights that JMJD6 confers an epigenetic vulnerability and mediates sunitinib sensitivity in renal cell carcinoma. <i>Clin. Transl. Med.</i> doi: 10.1002/ctm2.328.
RIME	Estrogen Receptor- $\alpha$	Mol. Cancer Res.	2021	Joseph L. Sottnik <i>et al.</i> Mediator of DNA Damage Checkpoint 1 (MDC1) Is a Novel Estrogen Receptor Coregulator in Invasive Lobular Carcinoma of the Breast. <i>Mol. Cancer Res.</i> doi: 10.1158/1541-7786.MCR-21-0025.
ChIP-Seq	Histone PTMs	Toxicol. Sci.	2021	Jo Jongpyo Lim <i>et al.</i> Neonatal Exposure to BPA, BDE-99, and PCB Produces Persistent Changes in Hepatic Transcriptome Associated With Gut Dysbiosis in Adult Mouse Livers. <i>Toxicol. Sci.</i> doi: 10.1093/toxsci/kfab104.
ChIP-qPCR	WT1	Molecular Oncology	2021	Alain Chebly <i>et al.</i> Exploring hTERT promoter methylation in cutaneous T-cell lymphomas. <i>Mol. Oncol.</i> doi: 10.1002/1878-0261.12946.
ATAC-Seq	NA	Blood	2021	Julie Agopian <i>et al.</i> GlcNAc is a mast-cell chromatin-remodeling oncometabolite that promotes systemic mastocytosis aggressiveness. <i>Blood.</i> doi: 10.1182/blood.2020008948.
ChIP-Seq	H3K27ac, RelA	Acta Neuropathologica	2021	Daniela Lotsch <i>et al.</i> Targeting fibroblast growth factor receptors to combat aggressive ependymoma. <i>Acta Neuropathol.</i> doi: 10.1007/s00401-021-02327-x.
ChIP-qPCR	H3K27ac	Molecular Cell	2021	Liang Ma <i>et al.</i> Co-condensation between transcription factor and coactivator p300 modulates transcriptional bursting kinetics. <i>Mol Cell.</i> doi: 10.1016/j.molcel.2021.01.031
ATAC-Seq, RNA-Seq	—	Cancer Immunol. Res.	2021	Justin C. Boucher <i>et al.</i> CD28 Costimulatory Domain-Targeted Mutations Enhance Chimeric Antigen Receptor T-cell Function. <i>Cancer Immunol. Res.</i> doi: 10.1158/2326-6066.CIR-20-0253.
RNA-Seq	—	Am. J. Physiol. Heart Circ. Physiol.	2021	Satvik Mareedu <i>et al.</i> Sarcolipin haploinsufficiency prevents dystrophic cardiomyopathy in mdx mice. <i>Am J Physiol Heart Circ Physiol.</i> doi: 10.1152/ajpheart.00601.2020.
ChIP-Seq	BRD4	Nature Chemical Biology	2021	Gabi Schutzius <i>et al.</i> BET bromodomain inhibitors regulate keratinocyte plasticity. <i>Nat. Chem. Biol.</i> doi: 10.1038/s41589-020-00716-z.
ChIP-qPCR	FOXO3	J. Cell Physiol.	2021	Siyun Wang <i>et al.</i> cMET promotes metastasis and epithelial-mesenchymal transition in colorectal carcinoma by repressing RKIP. <i>J. Cell Physiol.</i> doi: 10.1002/jcp.30142.
ATAC-Seq	—	Cancer Discovery	2021	Sonia Iyer <i>et al.</i> Genetically Defined Syngeneic Mouse Models of Ovarian Cancer as Tools for the Discovery of Combination Immunotherapy. <i>Cancer Discov.</i> doi: doi: 10.1158/2159-8290

Technique	Target	Journal	Year	Reference
ChIP-Seq	BRD2, BRD4, RNA Pol II	J. Immunol.	2021	Vinod Krishna <i>et al.</i> Integration of the Transcriptome and Genome-Wide Landscape of BRD2 and BRD4 Binding Motifs Identifies Key Superenhancer Genes and Reveals the Mechanism of Bet Inhibitor Action in Rheumatoid Arthritis Synovial Fibroblasts. <i>J. Immunol.</i> doi: doi: 10.4049/jimmunol.2000286.
ChIP-Seq	Multi-Histones	Cancer Cell	2021	Daniel Haag <i>et al.</i> . H3.3-K27M drives neural stem cell-specific gliomagenesis in a human iPSC-derived model. <i>Cancer Cell.</i> doi: 10.1016/j.ccell.2021.01.005.
ATAC-Seq	—	Cell Reports	2021	Ho Sun Jung <i>et al.</i> SOX17 integrates HOXA and arterial programs in hemogenic endothelium to drive definitive lympho-myeloid hematopoiesis. <i>Cell Rep.</i> doi: 10.1016/j.celrep.2021.108758.
ATAC-Seq	—	Cancers (Basel)	2021	Dongho Kim <i>et al.</i> Metformin Reduces Histone H3K4me3 at the Promoter Regions of Positive Cell Cycle Regulatory Genes in Lung Cancer Cells. <i>Cancers (Basel).</i> doi: 10.3390/cancers13040739.
ATAC-Seq	—	PLoS Pathog.	2021	Naveen Parmer <i>et al.</i> Intestinal-epithelial LSD1 controls goblet cell maturation and effector responses required for gut immunity to bacterial and helminth infection. <i>PLoS Pathog.</i> doi: 10.1371/journal.ppat.1009476.
PTM Quantitation	—	eLife	2021	Helen M. Tauc <i>et al.</i> Age-related changes in polycomb gene regulation disrupt lineage fidelity in intestinal stem cells. <i>eLife.</i> doi: 10.7554/eLife.62250
ChIP-Seq	H3K79me2	Diabetes	2021	Lin Shuai <i>et al.</i> DOT1L Regulates Thermogenic Adipocyte Differentiation and Function via Modulating H3K79 Methylation. <i>Diabetes.</i> doi: 10.2337/db20-1110
ChIP-Seq	TDF4	Cell Chem. Biol.	2021	Jiongjia Cheng <i>et al.</i> Small-molecule probe reveals a kinase cascade that links stress signaling to TCF/LEF and Wnt responsiveness. <i>Cell Chem. Biol.</i> doi: 10.1016/j.chembiol.2021.01.001
Mod Spec	—	Cancers (Basel)	2021	Anke Koeniger <i>et al.</i> Activation of Cilia-Independent Hedgehog/GLI1 Signaling as a Novel Concept for Neuroblastoma Therapy. <i>Cancers (Basel).</i> doi: 10.3390/cancers13081908
Histone PTM Analysis	—	Developmental Cell	2021	Igor L. Bado <i>et al.</i> The bone microenvironment increases phenotypic plasticity of ER + breast cancer cells. <i>Dev. Cell.</i> doi: 10.1016/j.devcel.2021.03.008
ATAC-Seq	—	Int. J. Mol. Sci.	2021	Lilas Courtot <i>et al.</i> Low Replicative Stress Triggers Cell-Type Specific Inheritable Advanced Replication Timing. <i>Int. J. Mol. Sci.</i> doi.org/10.3390/ijms22094959
RIME	RelA	Cell Reports	2021	James A. Heslop <i>et al.</i> GATA6 defines endoderm fate by controlling chromatin accessibility during differentiation of human-induced pluripotent stem cells. <i>Cell Rep.</i> doi: 10.1016/j.celrep.2021.109145
ChIP	H3K27ac, RelA	Acta Neuropathologica	2021	Daniela Lotsch <i>et al.</i> Targeting fibroblast growth factor receptors to combat aggressive ependymoma. <i>Acta Neuro.</i> doi: 10.1007/s00401-021-02327
ATAC-Seq	—	Nature Communications	2021	Joana Esteves de Lima <i>et al.</i> HIRA stabilizes skeletal muscle lineage identity. <i>Nature.</i> doi: 10.1038/s41467-021-23775-9

Technique	Target	Journal	Year	Reference
ChIP-Seq	ZNF263	ERJ Open Res.	2021	Sai Sneha Priya Nemani <i>et al.</i> COL4A3 expression in asthmatic epithelium depends on intronic methylation and ZNF263 binding. <i>ERJ Open Res.</i> doi: 10.1183/23120541.00802-2020
ChIP-Seq	BRD4, H3K27ac, H3K4me1	Mol. Ther. Methods Clin. Dev.	2021	Qingwei Wang <i>et al.</i> A hierarchical and collaborative BRD4/CEBPD partnership governs vascular smooth muscle cell inflammation. <i>Mol. Ther. Methods Clin. Dev.</i> doi: 10.1016/j.omtm.2021.02.021
ChIP-Seq	H2AK119Ub, Ring1b	Nature Chemical Biology	2021	Shirish Shukla <i>et al.</i> Small-molecule inhibitors targeting Polycomb repressive complex 1 RING domain. <i>Nature Chemical Biology.</i> doi: 10.1038/s41589-021-00815-5
ATAC-Seq	—	J. Mol. and Cell. Cardiology	2021	Douglas Chapski <i>et al.</i> Early adaptive chromatin remodeling events precede pathologic phenotypes and are reinforced in the failing heart. <i>J. Mol. and Cell. Cardiology.</i> doi: 10.1016/j.yjmcc.2021.07.002
RRBS	—	Nucleic Acids Research	2021	Miriam Recalde <i>et al.</i> The splicing regulator SLU7 is required to preserve DNMT1 protein stability and DNA methylation. <i>Nuc. Acids Res.</i> doi: 10.1093/narlgkab649
ATAC-Seq	—	EMBO Reports	2021	Raquel Sales-Gil <i>et al.</i> Non-redundant functions of H2A.Z.1 and H2A.Z.2 in chromosome segregation and cell cycle progression. <i>EMBO Reports.</i> doi: 10.15252/embr.202052061
RNA-Seq	—	Cells	2021	Melanie Lavaud <i>et al.</i> Overexpression of the Ubiquitin Specific Proteases USP43, USP41, USP27x and USP6 in Osteosarcoma Cell Lines: Inhibition of Osteosarcoma Tumor Growth and Lung Metastasis Development by the USP Antagonist PR619. <i>Cells.</i> doi: 10.33990/cells10092268
ChIP-Seq	FANCD2	Communications Biology	2021	Philippe Fernandes <i>et al.</i> FANCD2 modulates the mitochondrial stress response to prevent common fragile site instability. <i>Commun Biol.</i> doi: 10.1038/s42003-021-01647-8
RRBS	—	Science Advances	2021	Rina Baba <i>et al.</i> LSD1 enzyme inhibitor TAK-418 unlocks aberrant epigenetic machinery and improves autism symptoms in neurodevelopmental disorder models. <i>Sci Adv.</i> doi: 10.1126/sciadv.aba1187
RNA-Seq	—	Development	2021	Matthias Godart <i>et al.</i> The murine intestinal stem cells are highly sensitive to the modulation of the T3/TRa1-dependent pathway. <i>Development.</i> doi: 10.1242/dev.194357
ATAC-Seq	—	Nature Cancer	2021	Katherine R. Morel <i>et al.</i> EZH2 inhibition activates a dsRNA–STING–interferon stress axis that potentiates response to PD-1 checkpoint blockade in prostate cancer. <i>Nat Cancer.</i> doi: 10.1038/s43018-021-00185-w
ChIP-Seq	NR2F2	Sexual Development	2021	Fei Zhao <i>et al.</i> Molecular Actions Underlying Wolffian Duct Regression in Sexual Differentiation of Murine Reproductive Tracts. <i>Sex Dev.</i> doi: 10.1159/000513878
ChIP-Seq	H3K27ac	Mol. Neurobiol.	2020	Sunil Bhattarai <i>et al.</i> Modulation of Brain Pathology by Enhancer RNAs in Cerebral Ischemia. <i>Mol. Neurobiol.</i> doi: 10.1007/s12035-020-02194-9
ChIP-Seq	H3K9me3, H3K27me3	Arch. Toxicol.	2020	Masaki Fujioka <i>et al.</i> Dimethylarsinic acid (DMA) enhanced lung carcinogenesis via histone H3K9 modification in a transplacental mouse model. <i>Arch. Toxicol.</i> doi: 10.1007/s00204-020-02665-x.

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27me3	Cancer Cell	2020	Huairui Yuan <i>et al.</i> SETD2 Restricts Prostate Cancer Metastasis by Integrating EZH2 and AMPK Signaling Pathways. <i>Cancer Cell</i> . doi: 10.1016/j.ccell.2020.05.022
RIME	Bcl-6	Cell	2020	Leandro Venturutti <i>et al.</i> TBL1XR1 Mutations Drive Extranodal Lymphoma by Inducing a Pro-tumorigenic Memory Fate. <i>Cell</i> . doi: 10.1016/j.cell.2020.05.049.
RIME	Estrogen Receptor- $\alpha$	Sci Signal.	2020	Yin Li <i>et al.</i> A mutant form of ER $\alpha$ associated with estrogen insensitivity affects the coupling between ligand binding and coactivator recruitment. <i>Sci. Signal.</i> doi: 10.1126/scisignal.aaw4653.
ChIP-Seq	Progesterone Receptor	Sci. Signal.	2020	Margeaux Wetendorf <i>et al.</i> Constitutive expression of progesterone receptor isoforms promotes the development of hormone-dependent ovarian neoplasms. <i>Sci. Signal.</i> doi: 10.1126/scisignal.aaz9646.
ChIP-Seq	CEBP, H3K4me3, H3K27ac	Cell	2020	Ada Ndoja <i>et al.</i> Ubiquitin Ligase COP1 Suppresses Neuroinflammation by Degrading c/EBP $\beta$ in Microglia. <i>Cell</i> . doi: 10.1016/j.cell.2020.07.011.
ATAC-Seq	—	Nature Genetics	2020	Mark M. Pomerantz <i>et al.</i> Prostate cancer reactivates developmental epigenomic programs during metastatic progression. <i>Nat. Genet.</i> doi: 10.1038/s41588-020-0664-8.
ChIP-Seq	IRF-2	Nat. Cell Biol.	2020	Taku Sato <i>et al.</i> Regulated IFN signalling preserves the stemness of intestinal stem cells by restricting differentiation into secretory-cell lineages. <i>Nat. Cell Biol.</i> doi: 10.1038/s41556-020-0545-5.
ChIP-qPCR	BRD2, BRD4	Cardiovasc. Ther.	2020	Sylwia Wasiak <i>et al.</i> Epigenetic Modulation by Apabetalone Counters Cytokine-Driven Acute Phase Response In Vitro, in Mice and in Patients with Cardiovascular Disease. <i>Cardiovasc. Ther.</i> doi: 10.1155/2020/9397109.
ChIP-qPCR	BRG1, MITF	Mol. Cancer Ther.	2020	Florencia Rago <i>et al.</i> The Discovery of SWI/SNF Chromatin Remodeling Activity as a Novel and Targetable Dependency in Uveal Melanoma. <i>Mol. Cancer Ther.</i> doi: 10.1158/1535-7163
RNA-Seq	—	Nature	2020	Ana P. Gomes <i>et al.</i> Age-induced accumulation of methylmalonic acid promotes tumour progression. <i>Nature</i> . doi: 10.1038/s41586-020-2630-0.
ChIP-Seq	RNA Pol II	PLoS Genetics	2020	Katherine I. Farley-Barnes <i>et al.</i> Paired Box 9 (PAX9), the RNA polymerase II transcription factor, regulates human ribosome biogenesis and craniofacial development. <i>PLoS Genet.</i> doi: 10.1371/journal.pgen.1008967.
ChIP-Seq	H3K27ac	Transl. Oncol.	2020	Jie Cui <i>et al.</i> A zinc finger family protein, ZNF263, promotes hepatocellular carcinoma resistance to apoptosis via activation of ER stress-dependent autophagy. <i>Transl. Oncol.</i> doi: 10.1016/j.tranon.2020.100851.
RNA-Seq	—	Blood Adv.	2020	Iosifina P. Foskolou <i>et al.</i> The S enantiomer of 2-hydroxyglutarate increases central memory CD8 populations and improves CAR-T therapy outcome. <i>Blood Adv.</i> doi: 10.1182/bloodadvances.2020002309.
ChIP-Seq	b-catenin	Exp. Mol. Med.	2020	Jung-Yoon Yoo <i>et al.</i> $\beta$ -catenin activates TGF- $\beta$ -induced epithelial-mesenchymal transition in adenomyosis. <i>Exp. Mol. Med.</i> doi: 10.1038/s12276-020-00514-6
ChIP-Seq	Estrogen Receptor	FASEB J.	2020	Yukitomo Arao <i>et al.</i> The genomic regulatory elements for estrogen receptor alpha transactivation-function-1 regulated genes. <i>FASEB J.</i> doi: 10.1096/fj.202001435R.

Technique	Target	Journal	Year	Reference
ChIP-Seq	—	Nature Communications	2020	Gulfem D. Guler <i>et al.</i> Detection of early stage pancreatic cancer using 5-hydroxymethylcytosine signatures in circulating cell free DNA. <i>Nat. Commun.</i> doi: 10.1038/s41467-020-18965-w.
ChIP-Seq	SMARCA2, SMARCA4	Nature Communications	2020	Tharu M. Fernando <i>et al.</i> Functional characterization of SMARCA4 variants identified by targeted exome-sequencing of 131,668 cancer patients. <i>Nat. Commun.</i> doi: 10.1038/s41467-020-19402-8.
Mod Spec	—	Blood Adv.	2020	Tiziana Bruno <i>et al.</i> Che-1/AATF-induced transcriptionally active chromatin promotes cell proliferation in multiple myeloma. <i>Blood Adv.</i> doi: 10.1182/bloodadvances.2020002566.
RNA-Seq	—	Cancers (Basel)	2020	Mathilde Mullard <i>et al.</i> Sonic Hedgehog Signature in Pediatric Primary Bone Tumors: Effects of the GLI Antagonist GANT61 on Ewing's Sarcoma Tumor Growth. <i>Cancers (Basel)</i> . doi: 10.3390/cancers12113438.
RNA-Seq	—	iScience	2020	Aarti Mishra <i>et al.</i> Dynamic Neuroimmune Profile during Mid-life Aging in the Female Brain and Implications for Alzheimer Risk. <i>iScience</i> . doi: 10.1016/j.isci.2020.101829.
ChIP-Seq	H3K27ac, H3K27me3, H3K3me1, H3K9me3, H3K36me3	Nature Communications	2020	Yonghe Wu <i>et al.</i> Glioblastoma epigenome profiling identifies SOX10 as a master regulator of molecular tumour subtype. <i>Nat. Commun.</i> doi: 10.1038/s41467-020-20225-w.
RNA-Seq	—	Cancers (Basel)	2020	Sarah Morice <i>et al.</i> The YAP/TEAD Axis as a New Therapeutic Target in Osteosarcoma: Effect of Verteporfin and CA3 on Primary Tumor Growth. <i>Cancers (Basel)</i> . doi: 10.3390/cancers12123847.
ATAC-Seq	—	Nature	2020	Biraj Mahato <i>et al.</i> Pharmacologic fibroblast reprogramming into photoreceptors restores vision. <i>Nature</i> . 2020 doi: 10.1038/s41586-020-2201-4
ChIP-Seq	N-MYC, BRD4, H3K9me3, RNA Pol II	Nature Communications	2020	Maged Zeineldin <i>et al.</i> MYCN amplification and ATRX mutations are incompatible in neuroblastoma. <i>Nat. Commun.</i> doi: 10.1038/s41467-020-14682-6
ChIP-qPCR	HIF1alpha	Cell Reports	2020	Louise Ménégaut <i>et al.</i> Interplay between Liver X Receptor and Hypoxia Inducible Factor 1a Potentiates Interleukin-1b Production in Human Macrophages. <i>Cell Reports</i> . doi: 10.1016/j.celrep.2020.107665
ChIP-Seq	FLAG-tagged MRG15	Nature Metabolism	2020	Yuda Wei <i>et al.</i> MRG15 orchestrates rhythmic epigenomic remodelling and controls hepatic lipid metabolism. <i>Nat Metab.</i> 2020 doi: 10.1038/s42255-020-0203-z
Hi-C	—	Journal of Biological Chemistry	2020	Sylvia Hewitt <i>et al.</i> Estrogen receptor $\alpha$ (ER $\alpha$ )-binding super enhancers drive key mediators that control uterine estrogen responses in mice. <i>JBC</i> . 2020 doi: 10.1074/jbc.RA120.013666
ATAC-Seq	—	PNAS	2020	Kruttika Bhat <i>et al.</i> The dopamine receptor antagonist trifluoperazine prevents phenotype conversion and improves survival in mouse models of glioblastoma. <i>PNAS</i> . 2020 doi: 10.1073/pnas.1920154117



Technique	Target	Journal	Year	Reference
ATAC-Seq, ChIP-Seq	N-MYC, C-MYC, H3K4me1, H3K4me3, H3K27ac, H3K27me3	Scientific Data	2020	Kristen Upton <i>et al.</i> Epigenomic profiling of neuroblastoma cell lines. <i>Sci Data</i> . 2020 doi: 10.6084/m9.figshare.11993883
ChIP-Seq	H3K27ac, AFF4	Communications Biology	2020	Yujing Gao <i>et al.</i> Acetylation of histone H3K27 signals the transcriptional elongation for estrogen receptor alpha. <i>Commun Biol</i> . 2020 doi: 10.1038/s42003-020-0898-0
ChIP-Seq	HOXB13, H3K27ac, H3K27me3, H3K4me3, RNA Pol II	Experimental Cell Research	2020	Paul-Joseph Aspuria <i>et al.</i> HOXB13 controls cell state through super-enhancers. <i>Exp Cell Res</i> . 2020 doi: 10.1016/j.yexcr.2020.112039
ChIP-Seq	HA-tagged AHR1	mBio	2020	Sofia Ruben <i>et al.</i> Ahr1 and Tup1 Contribute to the Transcriptional Control of Virulence-Associated Genes in <i>Candida albicans</i> . <i>mBio</i> . 2020 doi: 10.1128/mBio.00206-20
ATAC-Seq	—	Cancer Discovery	2020	Di Zhao <i>et al.</i> Chromatin Regulator, CHD1, Remodels the Immunosuppressive Tumor Microenvironment in PTEN-Deficient Prostate Cancer. <i>Cancer Discov</i> . 2020 doi: 10.1158/2159-8290.CD-19-1352
ATAC-Seq	—	Diabetes	2020	Marion de Toledo <i>et al.</i> Lamin C counteracts glucose intolerance in aging, obesity and diabetes. <i>Diabetes</i> . 2020 doi: 10.2337/db19-0377.
ChIP-Seq	N-MYC, H3K27ac	Neoplasia	2020	Dongdong Chen <i>et al.</i> LIN28B promotes neuroblastoma metastasis and regulates PDZ binding kinase. <i>Neoplasia</i> . 2020 doi: 10.1016/j.neo.2020.04.001
ChIP-Seq	N-MYC, C-MYC	Frontiers in Oncology	2020	Robyn T. Sussman <i>et al.</i> CAMKV Is a Candidate Immunotherapeutic Target in MYCN Amplified Neuroblastoma. <i>Front. Oncol</i> . 2020 doi: 10.3389/fonc.2020.00302
ChIP-Seq	H3K4me3, H3K9me3, H3K27me3, H3R8me2a, H4K20me3	Theranostics	2020	Qian Zhang <i>et al.</i> Mdig promotes oncogenic gene expression through antagonizing repressive histone methylation markers. <i>Theranostics</i> 2020; 10(2):602-614.
ChIP-Seq	H3K4me1, H3K4me3, H3K9me3, H3K27me3, H3K27ac, H3K36me3, BRD4, MED1, MYC	Haematologica	2020	Aneta Mikulasova <i>et al.</i> Microhomology-mediated end joining drives complex rearrangements and over expression of MYC and PVT1 in multiple myeloma. <i>Haematologica</i> .doi:10.3324/haematol.2019.217927
ChIP-Seq	Androgen Receptor	Journal of Molecular Medicine	2019	Kinza Younas <i>et al.</i> Delayed endometrial decidualisation in polycystic ovary syndrome; the role of AR-MAGEA11. <i>J Mol Med</i> 2019 doi: 10.1007/s00109-019-01809-6
Mod Spec	—	Cell Reports	2019	Banushree Kumar <i>et al.</i> Quantitative Multiplexed ChIP Reveals Global Alterations that Shape Promoter Bivalency in Ground State Embryonic Stem Cells. <i>Cell Reports</i> . 28: 3274–3284.
RNA-Seq	—	The EMBO Journal	2019	Kathryn A. Jacobs <i>et al.</i> Paracaspase MALT1 regulates glioma cell survival by controlling endo-lysosome homeostasis. <i>EMBO J</i> . doi: 10.15252/embj.2019102030
ChIP-Seq	Estrogen Receptor	Cell	2019	Jane Guan <i>et al.</i> Therapeutic ligands antagonize estrogen receptor function by impairing its mobility. <i>Cell</i> . 178: 15

Technique	Target	Journal	Year	Reference
ATAC-Seq, ChIP-Seq	H3K4me3, H3K27ac, H3K27me3	Nature	2019	Nina Oberbeck <i>et al.</i> The RIPK4–IRF6 signalling axis safeguards epidermal differentiation and barrier function. <i>Nature</i> . 574: 249
ATAC-Seq	—	Cancer Cell	2019	Ana P. Gomes <i>et al.</i> Dynamic incorporation of histone H3 variants into chromatin is essential for acquisition of aggressive traits and metastatic colonization. <i>Cancer Cell</i> . 36: 402
ChIP-Seq	H3K27me3	Cancer Cell	2019	Zulekha A. Qadeer <i>et al.</i> ATRX in-frame fusion neuroblastoma is sensitive to EZH2 inhibition via modulation of neuronal gene signatures. <i>Cancer Cell</i> . 36: 512
ChIP-Seq	YAP1, NFIA	Nature Communications	2019	Kristian W. Pajtler <i>et al.</i> YAP1 subgroup supratentorial ependymoma requires TEAD and nuclear factor I-mediated transcriptional programmes for tumorigenesis. <i>Nat Commun</i> . 10: 3914
ChIP-Seq	BRG1, H3K9Ac	Nature Communications	2019	Min Liu <i>et al.</i> BRG1 attenuates colonic inflammation and tumorigenesis through autophagy-dependent oxidative stress sequestration. <i>Nat Commun</i> . 10: 4614
ChIP-Seq	BRD4	Nature Communications	2019	Nicolas Mercado <i>et al.</i> IRF2 is a master regulator of human keratinocyte stem cell fate. <i>Nat Commun</i> . 10: 4676
ChIP-Seq	RUNX1	Nature Communications	2019	Barbara Nicol <i>et al.</i> RUNX1 maintains the identity of the fetal ovary through an interplay with FOXL2. <i>Nat. Commun</i> . 10: 5116
ChIP-Seq	H3K27ac	Nature Communications	2019	Jayaram Vijayakrishnan <i>et al.</i> Identification of four novel associations for B-cell acute lymphoblastic leukaemia risk. <i>Nat Commun</i> . 10: 5348
ChIP-Seq	BRD4, H3K9me3	Neuron	2019	Jackie L. Norrie <i>et al.</i> Nucleome dynamics during retinal development. <i>Neuron</i> . 104:1
ChIP-qPCR	NRF2	Nature Immunology	2019	Anne-Valerie Burgener <i>et al.</i> SDHA gain-of-function engages inflammatory mitochondrial retrograde signaling via KEAP1–Nrf2. <i>Nat Immunol</i> . 20: 1311
ChIP-Seq, ChIP-qPCR	Androgen Receptor, Estrogen Receptor	Cell Reports	2019	Karyn Schmidt <i>et al.</i> The lncRNA SLNCR Recruits the Androgen Receptor to EGR1-Bound Genes in Melanoma and Inhibits Expression of Tumor Suppressor p21. <i>Cell Reports</i> . 27: 2493
ChIP-Seq	H4 panAc	Cell Metabolism	2019	Ramon C. Sun <i>et al.</i> Nuclear glycogenolysis modulates histone acetylation in human non-small cell lung cancers. <i>Cell Metab</i> . 30: 1
ChIP-Seq	RNA Pol II	Cell Stem Cells	2019	Kimberley N. Babos <i>et al.</i> Mitigating antagonism between transcription and proliferation allows near-deterministic cellular reprogramming. <i>Cell Stem Cell</i> . 25: 486
ChIP-Seq, RNA-Seq Data Analysis	ASH2L, H3K4me3	Cell Reports	2019	Liang Li <i>et al.</i> The COMPASS Family Protein ASH2L Mediates Corticogenesis via Transcriptional Regulation of Wnt Signaling. <i>Cell Reports</i> . 28: 698
Mod Spec	—	Nature Genetics	2019	Sara Martire <i>et al.</i> Phosphorylation of histone H3.3 at serine 31 promotes p300 activity and enhancer acetylation. <i>Nat. Genet</i> . 51: 941
ChIP-Seq	H3K27ac	Nature Communications	2019	Liqing Tian <i>et al.</i> Long-read sequencing unveils IGH-DUX4 translocation into the silenced IGH allele in B-cell acute lymphoblastic leukemia. <i>Nat. Commun</i> . 10: 2789

Technique	Target	Journal	Year	Reference
ChIP-Seq	SALL3, DNMT3B	Nature Communications	2019	Takuya Kuroda <i>et al.</i> SALL3 expression balance underlies lineage biases in human induced pluripotent stem cell differentiation. <i>Nat Commun.</i> 10: 2175
ChIP-Seq	H3K27me3, H3K27ac, H3K4me3	Investigative Ophthalmology and Visual Science	2019	Christina F. Tingle <i>et al.</i> Paradoxical Changes Underscore Epigenetic Reprogramming During Adult Zebrafish Extraocular Muscle Regeneration. <i>Invest Ophthalmol Vis Sci.</i> 2019;60:4991–4999
ChIP-Seq, ChIP-qPCR	Progesterone Receptor	The Journal of Clinical Endocrinology & Metabolism	2019	Ru-pin Alicia Chi <i>et al.</i> Human endometrial transcriptome and progesterone receptor cistrome reveal important pathways and epithelial regulators. <i>J Clin Endocrinol Metab.</i> doi: 10.1210/clinem/dgz117
ATAC-Seq	—	Haematologica	2019	James Ropa <i>et al.</i> SETDB1 mediated histone H3 lysine 9 methylation suppresses MLL-fusion target expression and leukemic transformation. <i>Haematologica.</i> doi: 10.3324/haematol.2019.223883
ChIP-Seq, Spike-in	H3K27me3	Haematologica	2019	Shao Xie <i>et al.</i> EZH2 inhibitors abrogate upregulation of trimethylation of H3K27 by CDK9 inhibitors and potentiate its activity against diffuse large B-cell lymphoma. <i>Haematologica.</i> doi:10.3324/haematol.2019.222935
ChIP-Seq, ChIP-qPCR	Androgen Receptor	Disease Models & Mechanisms	2019	Claire Nash <i>et al.</i> Genome-wide analysis of AR binding and comparison with transcript expression in primary human fetal prostate fibroblasts and cancer associated fibroblasts. <i>Mol Cell Endocrinol.</i> 471: 1
ChIP-Seq	H3K4me1, H3K4me3, H3K27ac, H3K27me3	BMC Biology	2019	Lingzhao Fang <i>et al.</i> Functional annotation of the cattle genome through systematic discovery and characterization of chromatin states and butyrate-induced variations. <i>BMC Biol.</i> 17: 68
ChIP-qPCR	BRD4, RelA	Clinical Epigenetics	2019	Laura Tsujikawa <i>et al.</i> Apabetalone (RVX-208) reduces vascular inflammation in vitro and in CVD patients by a BET-dependent epigenetic mechanism. <i>Clin Epigenet.</i> 11: 102
RIME	PPARb/d	Nucleic Acids Research	2019	Nathalie Legrand <i>et al.</i> PPARb/d recruits NCOR and regulates transcription reinitiation of ANGPTL4. <i>Nucleic Acids Res.</i> 47: 9573
ChIP-Seq	Progesterone Receptor, H3K27ac	Scientific Reports	2019	Diem T. Dinh <i>et al.</i> Tissue-specific progesterone receptor-chromatin binding and the regulation of progesterone-dependent gene expression. <i>Sci Rep.</i> 19: 9
Next-Gen Bisulfite-Seq	—	Plant Biotechnology Journal	2019	Aurine Verkest <i>et al.</i> Impact of differential DNA methylation on transgene expression in cotton ( <i>Gossypium hirsutum</i> L.) events generated by targeted sequence insertion. <i>Plant Biotechnol. J.</i> 17: 1236
ChIP-Seq	H3K9me3	Journal of Experimental Medicine	2019	Jun Hyung Sin <i>et al.</i> The epigenetic regulator ATF7ip inhibits IL2 expression, regulating Th17 responses. <i>J. Exp Med.</i> 216: 2024
ChIP-Seq	H3K27ac	Cancer Research	2019	Joesph Castillo <i>et al.</i> CBP/p300 drives the differentiation of Regulatory T cells through transcriptional and non-transcriptional mechanisms. <i>Cancer Res.</i> 79: 3916
ChIP-Seq, RNA-Seq	H3K4me3, H3K27me3	The Journal of Clinical Investigation	2019	Taiyi Huo <i>et al.</i> Induction of a cell-restricted Gc in dedifferentiating $\beta$ cells contributes to stress-induced $\beta$ -cell dysfunction. <i>J Clin Invest.</i> 5: 128351
ChIP-Seq	SMC1a, H3K4me1, H3K4me3, H3K27ac	Journal of Biological Chemistry	2019	Sylvia C. Hewitt <i>et al.</i> A distal super enhancer mediates estrogen-dependent mouse uterine-specific gene transcription of Igf1 (Insulin-like growth factor 1). <i>J Biol Chem.</i> 294: 9746

Technique	Target	Journal	Year	Reference
ChIP-Seq and ChIP-qPCR	CBP, p300, H3K27ac	Cell Reports	2019	Denise de Almeida Nagata <i>et al.</i> Regulation of Tumor-Associated Myeloid Cell Activity by CBP/EP300 Bromodomain Modulation of H3K27 Acetylation. <i>Cell Reports</i> . 27:269
ChIP-Seq, RNA-Seq	GR	Journal of the American Heart Association	2019	Elena Severinova <i>et al.</i> Glucocorticoid Receptor Binding and Transcriptome Signature in Cardiomyocytes. <i>J Am Heart Assoc</i> . 8:e011484. DOI: 10.1161/JAHA.118.011484.
ChIP-Seq	H4K20me1	Journal of Autoimmunity	2019	Keqi Fan <i>et al.</i> CRL4DCAF2 is required for mature T-cell expansion via Aurora B-regulated proteasome activity. <i>Journal of Autoimmunity</i> . 96:74
MeDIP-Seq	—	Epigenomes	2019	Liliana Ferreira <i>et al.</i> Uncovering Differentially Methylated Regions (DMRs) in a Salt-Tolerant Rice Variety under Stress: One Step towards New Regulatory Regions for Enhanced Salt Tolerance. <i>Epigenomes</i> . 3:4
ChIP-qPCR	ChREBP	Nature Metabolism	2019	Pauline Morigny <i>et al.</i> Interaction between hormone-sensitive lipase and ChREBP in fat cells controls insulin sensitivity. <i>Nature Metabolism</i> . 1: 133
ChIP-Seq	H3K4me1, H3K4me3, H3K27ac	Cell Metabolism	2019	Liming Du <i>et al.</i> IGF-2 Preprograms Maturing Macrophages to Acquire Oxidative Phosphorylation-Dependent Antiinflammatory Properties. <i>Cell Metabolism</i> . 29: 1
ChIP-Seq	RNA Pol II, H3K27ac	Journal of Neuro-Oncology	2019	Isabel Tegeder <i>et al.</i> Functional relevance of genes predicted to be affected by epigenetic alterations in atypical teratoid/rhabdoid tumors. <i>J. Neurooncol.</i> <a href="https://doi.org/10.1007/s11060-018-03018-6">https://doi.org/10.1007/s11060-018-03018-6</a>
ChIP-Seq	EGR1, H3K27ac	Journal of Steroid Biochemistry & Molecular Biology	2019	Maria Szwarc <i>et al.</i> Early growth response 1 transcriptionally primes the human endometrial stromal cell for decidualization. <i>J. Steroid Biochem. Mol. Biol.</i> <a href="https://doi.org/10.1016/j.jsbmb.2019.01.021">https://doi.org/10.1016/j.jsbmb.2019.01.021</a>
ChIP-Seq	H3K27me3, H3K4me3	Acta Neuropathologica	2019	Andre Silveira <i>et al.</i> H3.3 K27M depletion increases differentiation and extends latency of diffuse intrinsic pontine glioma growth in vivo. <i>Acta Neuropathol.</i> <a href="https://doi.org/10.1007/s00401-019-01975-4">https://doi.org/10.1007/s00401-019-01975-4</a>
ChIP-Seq	RNA Pol II	Agronomy	2019	Sonja Klemme <i>et al.</i> Selection of Salicylic Acid Tolerant Epilines in <i>Brassica napus</i> . <i>Agronomy</i> . 9: 92
ChIP-Seq	LSD1, H3K4me2, H3K27ac	Science Signaling	2019	Arnaud Augert <i>et al.</i> Targeting NOTCH activation in small cell lung cancer through LSD1 inhibition. <i>Sci. Signal.</i> 12, eaau2922
ChIP-qPCR	SRF and CEBP	Nature Communications	2019	Shilpita Sarcar <i>et al.</i> Next-generation muscle-directed gene therapy by in silico vector design. <i>Nat Commun.</i> <a href="https://doi.org/10.1038/s41467-018-08283-7">https://doi.org/10.1038/s41467-018-08283-7</a>
ChIP-Seq	H3K27ac, H3K4me3, CTCF, and NR4A3	Nature Communications	2019	Florian Haller <i>et al.</i> Enhancer hijacking activates oncogenic transcription factor NR4A3 in acinic cell carcinomas of the salivary glands. <i>Nat Commun.</i> 10: 336
ChIP-Seq	Gfi1, Lsd1	Nature Communications	2019	Catherine Lee <i>et al.</i> Lsd1 as a therapeutic target in Gfi1-activated medulloblastoma. <i>Nat Commun.</i> 10: 332

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27me3, H3K4me1, H3K4me3, H3K9me3, H3K36me3, EZH2, SMARCB1, SUZ12, SMARCA4, REST	Cancer Cell	2019	Serap Erkek <i>et al.</i> Comprehensive Analysis of Chromatin States in Atypical Teratoid/Rhabdoid Tumors Identified Diverging Roles for SWI/SNF and Polycomb in Gene Regulation. <i>Cancer Cell</i> . 35: 95
ChIP-Seq	BRG1	The Journal of Clinical Investigation	2019	Yufeng Ding <i>et al.</i> Chromatin remodeling ATPase BRG1 and PTEN are synthetic lethal in prostate cancer. <i>J Clin Invest</i> . <a href="https://doi.org/10.1172/JCI123557">https://doi.org/10.1172/JCI123557</a>
ChIP-qPCR	PPAR gamma	International Journal of Obesity	2019	Laura Butruille <i>et al.</i> Maternal high-fat diet during suckling programs visceral adiposity and epigenetic regulation of adipose tissue stearoyl-CoA desaturase-1 in offspring. <i>Int. J. Obes</i> . <a href="https://doi.org/10.1038/s41366-018-0310-z">https://doi.org/10.1038/s41366-018-0310-z</a>
ChIP-Seq	HA-tagged JP2NT, TBP, MEF2C	Science	2018	Ang Guo <i>et al.</i> E-C coupling structural protein junctophilin-2 encodes a stress-adaptive transcription regulator. <i>Science</i> . 362: 6421
ChIP-qPCR	p65/RelA	Human Molecular Genetics	2018	Jeffrey R. Gehlhausen <i>et al.</i> A proteasome-resistant fragment of NIK mediates oncogenic NF- $\kappa$ B signaling in schwannomas. <i>Clin Cancer Res</i> . 25:1601. doi: 10.1093/hmg/ddy361
ChIP-Seq	Nrf2	Journal of Biological Chemistry	2018	Junsheng Fu <i>et al.</i> Hyperactivity of the transcription factor Nrf2 causes metabolic reprogramming in mouse esophagus. <i>J. Biol. Chem</i> . 294: 327
ChIP-Seq	H3K36me2	Cell reports	2018	Jingjing Chen <i>et al.</i> Methyltransferase Nsd2 Ensures Germinal Center Selection by Promoting Adhesive Interactions between B Cells and Follicular Dendritic Cells. <i>Cell Reports</i> . 25: 3393
ChIP-Seq	H3K27ac	Clinical Cancer Research	2018	Shori Saito <i>et al.</i> Eradication of central nervous system leukemia of T-cell origin with a brain-permeable LSD1 inhibitor. <i>Clin Cancer Res</i> . 25:1601
ChIP-Seq	RNA Pol II, Brd4, H3K9me3	Cell Reports	2018	Lu Want <i>et al.</i> Retinal Cell Type DNA Methylation and Histone Modifications Predict Reprogramming Efficiency and Retinogenesis in 3D Organoid Cultures. <i>Cell Reports</i> . 22: 2601
ATAC-Seq	—	Cell Reports	2018	Xiaolong Zhang <i>et al.</i> OX40 Costimulation inhibits Foxp3 expression and Treg induction via BATF3-dependent and independent mechanisms. <i>Cell Reports</i> . 24: 607.
Mod Spec	—	Acta Neuropathologica	2018	Leah Katz <i>et al.</i> Loss of histone H3K27me3 identifies a subset of meningiomas with increased risk of recurrence. <i>Acta Neuropathol</i> . 135: 6.
RIME	TRPS1	Cell Reports	2018	Robert M. Witwicky <i>et al.</i> TRPS1 is a lineage-specific transcriptional dependency in breast cancer. <i>Cell</i> . 175: 1.
ChIP-Seq	H3K27ac	Cell	2018	Xun Huang <i>et al.</i> Targeting epigenetic crosstalk as a therapeutic strategy for EZH2-aberrant solid tumors. <i>Cell</i> . 175: 1.
ChIP-Seq	BRD4, RNA Pol II	Cancer Cell	2018	Elizabeth Stewart <i>et al.</i> Identification of therapeutic targets in rhabdomyosarcoma through integrated genomic, epigenomic, and proteomic analyses. <i>Cancer Cell</i> . 34: 1.
ChIP-Seq	Sox17	Nature Communications	2018	Xiaoqiu Wang <i>et al.</i> SOX17 regulates uterine epithelial-stromal cross-talk acting via a distal enhancer upstream of <i>Ihh</i> . <i>Nat Commun</i> . 9: 4421

Technique	Target	Journal	Year	Reference
ChIP-Seq, RNA-Seq	NRF2, RNA Pol II	Nature Communications	2018	David Olganier <i>et al.</i> Nrf2 negatively regulates STING indicating a link between antiviral sensing and metabolic reprogramming. <i>Nat Commun.</i> 9:3506
ChIP-Seq, RIME	p300, CBP, CDK9, BRD4, H3K27ac, H3K4me1, H3K4me3, H3K18Ac	Cell Reports	2018	Ryan Raisner <i>et al.</i> Enhancer activity requires CBP/P300 bromodomain-dependent histone H3K27 acetylation. <i>Cell Reports.</i> 24: 1722.
ChIP-Seq, RNA-Seq	H3K27ac, Med1, RNA Pol II	The Journal of Clinical Investigation	2018	Maria Florencia Martinez <i>et al.</i> Super-Enhancers maintain renin-expressing cell identity and memory to preserve multi-system homeostasis. <i>J. Clin Invest.</i> Advance Online Publication doi: 10.1172/JCI121361.
ChIP-Seq,	BRD4, H3K27ac	JCI Insight	2018	Anisley Valenciana <i>et al.</i> Transcriptional targeting of oncogene addiction in medullary thyroid cancer. <i>JCI Insight.</i> 3: e122225.
RNA-Seq	—	PAIN	2018	Pradipta Ray <i>et al.</i> Comparative transcriptome profiling of the human and mouse dorsal root ganglia: an RNA-Seq based resource for pain and sensory neuroscience research. <i>Pain.</i> 159: 7.
ChIP-Seq	ELL2	The Journal Of Immunology	2018	Ashley M. Nelson <i>et al.</i> RNA splicing in the transition from B cells to antibody-secreting cells: The influences of ELL2, small nuclear RNA, and endoplasmic reticulum stress. <i>J Immunol.</i> Advance online publication doi: 10.4049/jimmunol
ChIP-Seq	MTA1	Cancer Medicine	2018	Nasir A. Butt <i>et al.</i> Targeting MTA1/HIF-1 $\alpha$ signaling by pterostilbene in combination with histone deacetylase inhibitor attenuates prostate cancer progression. <i>Cancer Med.</i> 6: 2673
ChIP-Seq, Next-Gen Bisulfite-Seq	H3K4me3, RNA Pol II	Agronomy	2018	Martin Schmidt <i>et al.</i> Methylome and epialleles in rice epilines selected for energy use efficiency. <i>Agronomy.</i> 8: 163
ChIP-Seq	GATA2	The American Journal of Human Genetics	2018	Katelyn M. Mika <i>et al.</i> An ancient fecundability-associated polymorphism creates a GATA2 binding site in a distal enhancer of HLA-F. <i>Am J Hum Genet.</i> 103: 509
ChIP-Seq	FOXL2	Human Molecular Genetics	2018	Barbara Nicol <i>et al.</i> Genome-wide identification of FOXL2 binding and characterization of FOXL2 feminizing action in the fetal gonads. <i>Hum Mol Genet.</i> Advance online publication doi: 10.1093/hmg/ddy312
MeDIP-Seq	5-Methylcytosine	Scientific Reports	2018	Jose P. Silva <i>et al.</i> Analysis of diet-induced differential methylation, expression, and interactions of lncRNA and protein-coding genes in mouse liver. <i>Sci Rep.</i> 8: 11537
ChIP-Seq	ZFP24	Cell Reports	2018	Benayahu Elbaz <i>et al.</i> Phosphorylation state of ZFP24 controls oligodendrocyte differentiation. <i>Cell Reports.</i> 23: 2254
ChIP-Seq, Histone PTM	H3K9me3, H3K27me3	Cell Reports	2018	Jessica Camacho <i>et al.</i> The memory of environmental chemical exposure in <i>C. elegans</i> is dependent on the jumonji demethylases <i>jmjd-2</i> and <i>jmjd-3/utx-1</i> . <i>Cell Reports.</i> 23: 2392
ChIP-Seq	TET1, DNMT1	Nature Genetics	2018	Nipun Verma <i>et al.</i> TET proteins safeguard bivalent promoters from de novo methylation in human embryonic stem cells. <i>Nat Genet.</i> 50: 83
ChIP-Seq	EZH2	Journal of Experimental Medicine	2018	Xingli Zhang <i>et al.</i> Macrophage/microglial Ezh2 facilitates autoimmune inflammation through inhibition of Socs3. <i>J Exp Med.</i> 215: 1365

Technique	Target	Journal	Year	Reference
ChIP-Seq	STAT3,STAT5, ROCK2	Scientific Reports	2018	Wei Chen <i>et al.</i> ROCK2, but not ROCK1 interacts with phosphorylated STAT3 and co-occupies TH17/TFH gene promoters in TH17-activated human T cells. <i>Sci Rep.</i> 8: 16636
ChIP-Seq	Bhlhe40	PLoS One	2018	Kelly A. Hamilton <i>et al.</i> Mice lacking the transcriptional regulator Bhlhe40 have enhanced neuronal excitability and impaired synaptic plasticity in the hippocampus. <i>PLoS One.</i> 13: e0196223
ChIP-Seq	RNA Pol II	PLoS One	2018	Liana Basova <i>et al.</i> Dopamine and its receptors play a role in the modulation of CCR5 expression in innate immune cells following exposure to Methamphetamine: Implications to HIV infection. <i>PLoS One.</i> 13: e0199861
ChIP-qPCR	Glucocorticoid Receptor, TIF1	Arthritis & Rheumatology	2018	Yanhua Hu <i>et al.</i> Development of a molecular signature to monitor pharmacodynamic response mediated by in vivo administration of glucocorticoids. <i>Arthritis Rheumatol.</i> 70: 1331
ChIP-Seq	H3K27me3	Acta Neuropathologica Communications	2018	David Castel <i>et al.</i> Transcriptomic and epigenetic profiling of 'diffuse midline gliomas, H3 K27M-mutant' discriminate two subgroups based on the type of histone H3 mutated and not supratentorial or infratentorial location. <i>Acta Neuropathol Commun.</i> 6: 17
ChIP-Seq	NRF2	Molecular Pharmacology	2018	Rance Nault <i>et al.</i> Comparison of hepatic NRF2 and AHR binding in 2,3,7,8- tetrachlorodibenzo-p-dioxin (TCDD) treated mice demonstrates NRF2- independent PKM2 induction. <i>Mol Pharmacol.</i> 94: 876
ChIP-Seq	HDAC2, H3K27ac, H3K27me3	Cancer Immunology, Immunotherapy	2018	David Briere <i>et al.</i> The class I/IV HDAC inhibitor mocetinostat increases tumor antigen presentation, decreases immune suppressive cell types and augments checkpoint inhibitor therapy. <i>Cancer Immunol Immunother.</i> 67: 381
ChIP-Seq, ChIP-qPCR	TRRAP	Journal of Cell Biology	2018	Zhao Wang <i>et al.</i> TRRAP is a central regulator of human multiciliated cell formation. <i>J Cell Biol.</i> 217: 1941
ChIP-Seq	BRD4	Blood	2018	Rebecca Kohnken <i>et al.</i> Diminished microRNA-29b level is associated with BRD4 mediated activation of oncogenes in cutaneous T-cell lymphoma. <i>Blood.</i> 131: 771
RNA-Seq	—	Cell Communication and Signaling	2018	Keri Callegari <i>et al.</i> Pharmacological inhibition of LSD1 activity blocks REST-dependent medulloblastoma cell migration. <i>Cell Commun Signal.</i> 16: 60
ChIP-Seq	MTA1	Molecular Oncology	2018	Avinash Kumar <i>et al.</i> MTA drives malignant progression and bone metastasis in prostate cancer
ChIP-Seq	RNA Pol II, H2A.Z, CDK9, ANP32e	Biochimica et Biophysica Acta	2018	Hyewon Shin <i>et al.</i> Transcriptional regulation mediated by H2A.Z via ANP32e-dependent inhibition of protein phosphatase 2A. <i>Biochim Biophys Acta.</i> 1861: 481
ChIP-Seq	H3K4me3, H3K9me3, H3K27me3	Proceedings of the Royal Society B	2018	Theresa K. Kelly <i>et al.</i> Epigenetic regulation of transcriptional plasticity associated with developmental song learning. <i>Proc Biol Sci.</i> 285: 20180160
ChIP-Seq	SIRT1	Journal of Neuroimmune Pharmacology	2018	Nikki Bortell <i>et al.</i> Sirtuin 1-chromatin-binding dynamics points to a common mechanism regulating inflammatory targets in SIV infection and in the aging brain. <i>J Neuroimmune Pharmacol.</i> 13: 163

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27me3	G3 - Genes, Genomics Genetics	2018	James Ferguson <i>et al.</i> PRC2 is dispensable in vivo for b-catenin-mediated repression of chondrogenesis in the mouse embryonic cranial mesenchyme. <i>G3</i> . 8: 491
MeDIP-Seq	5-Methylcytosine	Nature	2017	Rama S. Akondy <i>et al.</i> Origin and differentiation of human memory CD8 T cells after vaccination. <i>Nature</i> . 552: 362
ChIP-Seq	10 different Histone mods plus CTCF, RNA Pol II, & BRD4	Nature	2017	Elizabeth Stewart <i>et al.</i> Orthotopic patient-derived xenografts of paediatric solid tumours. <i>Nature</i> . 549: 96
ChIP-Seq	BACH2	Nature Communications	2017	Nicolas Hipp <i>et al.</i> IL-2 imprints human naive B cell fate towards plasma cell through ERK/ELK1-mediated BACH2 repression. <i>Nat Commun</i> . 8: 1443
ChIP-qPCR	H3K9Ac, RNA Pol II, ChREBP, PPARa	Cell Reports	2017	Alison Iroz <i>et al.</i> A specific ChREBP and PPARa cross-talk is required for the glucose-mediated FGF21 response. <i>Cell Reports</i> . 21: 403
ChIP-Seq	H3K79me2	Molecular Cancer Therapy	2017	Carly T. Campbell <i>et al.</i> Mechanisms of Pinometostat (EPZ-5676) treatment-emergent resistance in MLL-rearranged leukemia. <i>Mol Cancer Ther</i> . 16: 1669
Next-Gen Bisulfite-Seq, ChIP-Seq	KLF4	Leukemia	2017	Y Shen <i>et al.</i> Inactivation of KLF4 promotes T-cell acute lymphoblastic leukemia and activates the MAP2K7 pathway. <i>Leukemia</i> . 31: 1314
ChIP-Seq	H3K27ac, H3K4me1	Epigenomics	2017	John P. Thomson <i>et al.</i> Defining baseline epigenetic landscapes in the rat liver. <i>Epigenomics</i> . 9: 1503
ChIP-Seq	H3K36me3	The Journal of Clinical Investigation	2017	Huairui Yuan <i>et al.</i> Histone methyltransferase SETD2 modulates alternative splicing to inhibit intestinal tumorigenesis. <i>J Clin Invest</i> . 127: 3375
RIME	Estrogen Receptor, Progesterone Receptor	Cancer Research	2017	Jessica Finlay-Schultz <i>et al.</i> Breast cancer suppression by progesterone receptors is mediated by their modulation of estrogen receptors and RNA polymerase III. <i>Cancer Res</i> . 77: 4934
ChIP-Seq	EZH2, H3K27me3, H3K4me3	Molecular Cancer Therapy	2017	Dorothy Brach <i>et al.</i> EZH2 inhibition by tazemetostat results in altered dependency on B-cell activation signaling in DLBCL. <i>Mol Cancer Ther</i> . 16:2586
ChIP-Seq Spike-in	p300, H3K27ac, Androgen Receptor	Cancer Research	2017	Lingyan Jin <i>et al.</i> Therapeutic targeting of the CBP/p300 bromodomain blocks the growth of castration-resistant prostate cancer. <i>Cancer Res</i> . 77: 5564
ChIP-qPCR	IRF5, NFkB (p65)	Journal of Biological Chemistry	2017	Leah Cushing <i>et al.</i> IRAK4 kinase activity controls Toll-like receptor induced inflammation through the transcription factor IRF5 in primary human monocytes. <i>J Biol Chem</i> . 292: 18689.
ChIP-Seq	RNA Pol II	Journal of Biological Chemistry	2017	Leah A. Gates <i>et al.</i> Acetylation on histone H3 lysine 9 mediates a switch from transcription initiation to elongation. <i>J Biol Chem</i> . 292: 14456
ChIP-Seq ChIP-qPCR	Estrogen Receptor	Endocrinology	2017	Sylvia C. Hewitt <i>et al.</i> Role of ERa in mediating female uterine transcriptional Responses to IGF1. <i>Endocrinology</i> . 158: 2427
ATAC-Seq ChIP-Seq	H3K4me3, H3K9me3	Cancer Cell	2017	Gulfem Dilek Guler <i>et al.</i> Repression of stress-induced LINE-1 expression protects cancer cell subpopulations from lethal drug exposure. <i>Cancer Cell</i> . 32: 1



Technique	Target	Journal	Year	Reference
ChIP-Seq	BRD4, H3K9me3, RNA Pol II	Neuron	2017	Issam Aldiri <i>et al.</i> The dynamic epigenetic landscape of the retina during development, reprogramming and tumorigenesis. <i>Neuron</i> . 94: 550
ChIP-Seq	Lmx1b	Development	2017	Endika Haro <i>et al.</i> Lmx1b-targeted cis-regulatory modules involved in limb dorsalization. <i>Dev</i> . Advance online publication doi: 10.1242/dev.146332
ChIP-Seq	EZH2	Proceedings of the Natl. Academy of Sciences	2017	Yongfeng Liu <i>et al.</i> Epithelial EZH2 serves as an epigenetic determinant in experimental colitis by inhibiting TNF $\alpha$ -mediated inflammation and apoptosis. <i>Proc Natl Acad Sci</i> . 114: E3796
ChIP-Seq	Estrogen Receptor	Endocrinology	2017	Laurel A. Coons <i>et al.</i> DNA sequence constraints define functionally active steroid nuclear receptor binding sites in chromatin. <i>Endocrinology</i> . Advance online publication doi: 10.1210/en.2017-00468
ChIP-Seq	ETV5	Proceedings of the Natl. Academy of Sciences	2017	Zhen Zhang <i>et al.</i> Transcription factor Etv5 is essential for the maintenance of alveolar type II cells. <i>Proc Natl Acad Sci</i> . 114: E3903
ChIP-Seq	Androgen Receptor	Molecular and Cellular Endocrinology	2017	Claire Nash <i>et al.</i> Genome-wide analysis of AR binding and comparison with transcript expression in primary human fetal prostate fibroblasts and cancer associated fibroblasts. <i>Mol Cell Endocrinol</i> . Advance online publication doi: 10.1016/j.mce.2017.05.006
ChIP-Seq Analysis	H3K27me3	Scientific Reports	2017	Brid O'Leary <i>et al.</i> Long non-coding RNA PARTICLE bridges histone and DNA methylation. <i>Sci Rep</i> . 7: 1790
ChIP-Seq	H3K36me2	The Journal of Clinical Investigation	2017	Ni Li <i>et al.</i> AKT-mediated stabilization of histone methyltransferase WHSC1 promotes prostate cancer metastasis. <i>J Clin Invest</i> . 127: 1284
ChIP-Seq ChIP-qPCR	EZH2, H3K27me3, H3K27ac	Cancer Cell	2017	Eric E. Gardner <i>et al.</i> Chemosensitive relapse in small cell lung cancer proceeds through an EZH2-SLFN11 axis. <i>Cancer Cell</i> . 31: 286
ChIP-Seq	TRIM28	Cell Reports	2017	Per Ludvik Brattas <i>et al.</i> TRIM28 controls a gene regulatory network based on endogenous retroviruses in human neural progenitor cells. <i>Cell Reports</i> 18: 1
ChIP-Seq	Cbx3/HP1	Scientific Reports	2017	Michael Sun <i>et al.</i> Cbx3/HP1g deficiency confers enhanced tumor-killing capacity on CD8 <sup>+</sup> T cells. <i>Sci Rep</i> . 7: 42888
ChIP-Seq	H3K27ac	Scientific Reports	2017	Yao Shen <i>et al.</i> Epigenetic and genetic dissections of UV-induced global gene dysregulation in skin cells through multi-omics analyses. <i>Sci Rep</i> . 7: 42646
ChIP-qPCR	SIRT1	Scientific Reports	2017	Jung-Yoon Yoo <i>et al.</i> KRAS activation and over-expression of SIRT1/BCL6 contributes to the pathogenesis of endometriosis and progesterone resistance. <i>Sci Rep</i> . 7: 6765
MeDIP-Seq	5-methylcytosine	Cerebral Cortex	2017	Daniela Grassi <i>et al.</i> Neuronal activity, TGF-Signaling and unpredictable chronic stress modulate transcription of Gadd45 family members and DNA methylation in the hippocampus. <i>Cereb Cortex</i> . 27: 4166
ChIP-Seq	SRC-2	PLoS Genetics	2017	Shruthy Suresh <i>et al.</i> SRC-2-mediated coactivation of anti-tumorigenic target genes suppresses MYC-induced liver cancer. <i>PLoS Genet</i> . 13: e1006650

Technique	Target	Journal	Year	Reference
ChIP-Seq	SLY1	Cell Death and Differentiation	2017	Charlotte Moretti <i>et al.</i> SLY regulates genes involved in chromatin remodeling and interacts with TBL1XR1 during sperm differentiation. <i>Cell Death Differ.</i> Advance online publication doi: 10.1038/cdd.2017.32
ChIP-Seq	G34R	Acta Neuropathologica Communications	2017	Farhana Haque <i>et al.</i> Evaluation of a novel antibody to define histone 3.3 G34R mutant brain tumours. <i>Acta Neuropathol Commun.</i> 5: 45
ChIP-Seq	SALL4	Journal of Hematology & Oncology	2017	Lina Yang <i>et al.</i> The stem cell factor SALL4 is an essential transcriptional regulator in mixed lineage leukemia-rearranged leukemogenesis. <i>J Hematol Oncol.</i> 10: 159
ChIP-qPCR	EZH2	Mechanisms of Ageing and Development	2017	Khyobeni Mozhui <i>et al.</i> Conserved effect of aging on DNA methylation and association with EZH2 polycomb protein in mice and humans. <i>Mech Ageing Dev.</i> 162: 27
ChIP-Seq Spike-in	H3K27me3, H3K4me3, H3K9me3	PLoS One	2016	Brian Egan <i>et al.</i> An alternative approach to ChIP-Seq normalization enables detection of genome-wide changes in histone H3 lysine 27 trimethylation upon EZH2 inhibition. <i>PLoS One.</i> 11: e0166438
Low Cell ChIP-Seq	H3K27me3, CTCF	Nature Genetics	2016	Joachim Weischenfeldt <i>et al.</i> Pan-cancer analysis of somatic copy-number alterations implicates IRS4 and IGF2 in enhancer hijacking. <i>Nat Genet.</i> Advance online publication doi: 10.1038/ng.3722
ChIP-Seq	OTX2, MITF, BRD4, H3K27ac	Cancer Cell	2016	Pascal D. Johann <i>et al.</i> Atypical teratoid/rhabdoid tumors are comprised of three epigenetic subgroups with distinct enhancer landscapes. <i>Cancer Cell.</i> 29: 379
ChIP-Seq	GATA2, Progesterone Receptor	Cell Reports	2016	Cory A. Rubel <i>et al.</i> A Gata2-dependent transcription network regulates uterine progesterone responsiveness and endometrial function. <i>Cell Reports.</i> 17: 1414
ChIP-Seq	Androgen Receptor	Molecular Cell	2016	Boyu Zhang <i>et al.</i> Non-Cell-Autonomous Regulation of Prostate Epithelial Homeostasis by Androgen Receptor. <i>Mol Cell.</i> 63: 976
ChIP-Seq Spike-in	BRD4	The Journal of Clinical Investigation	2016	Mark L. McClelland <i>et al.</i> CCAT1 is an enhancer-templated RNA that predicts BET sensitivity in colorectal cancer. <i>J Clin Invest.</i> 126: 639
hMeDIP-Seq		Nature Communications	2016	Carolina M. Greco <i>et al.</i> DNA hydroxymethylation controls cardiomyocyte gene expression in development and hypertrophy. <i>Nat Commun.</i> 7: 12418
ChIP-qPCR	Androgen Receptor	Cell Reports	2016	Karyn Schmidt <i>et al.</i> The lncRNA SLNCR1 mediates melanoma invasion through a conserved SRA1-like region. <i>Cell Rep.</i> 15: 2025
Low Cell ChIP-Seq	H3K4me3, H3K27ac, H3K27me3	Biology of Reproduction	2016	Kazadi Mutoji <i>et al.</i> TSPAN8 expression distinguishes spermatogonial stem cells in the prepubertal mouse testis. <i>Biol Reprod.</i> 95: 17
ChIP-Seq	HDAC1, H3K27ac	Cancer Discovery	2016	Anjali Mishra <i>et al.</i> Mechanism, consequences, and therapeutic targeting of abnormal IL15 signaling in cutaneous T-cell lymphoma. <i>Cancer Discov.</i> 6: 986
ChIP-Seq	V5-tagged DUX4	Human Molecular Genetics	2016	Jocelyn O. Eidahl <i>et al.</i> Mouse Dux is myotoxic and shares partial functional homology with its human paralog DUX4. <i>Hum Mol Genet.</i> Advance online publication doi: 10.1093/hmg/ddw287
ChIP-Seq	H3K27me3	Molecular and Cellular Biology	2016	Kyung Hyun Yoo <i>et al.</i> Histone demethylase KDM6A controls the mammary luminal lineage through enzyme-independent mechanisms. <i>Mol Cell Biol.</i> 36: 2108

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27ac	Nature Genetics	2016	Pratiti Bandopadhyay <i>et al.</i> MYB-QKI rearrangements in angiocentric glioma drive tumorigenicity through a tripartite mechanism. <i>Nat Genet.</i> 48: 273
ChIP-Seq Spike-in	KDM5C	Cell Reports	2016	Shigeki Iwase <i>et al.</i> A mouse model of X-linked intellectual disability associated with impaired removal of histone methylation. <i>Cell Reports.</i> 14: 1
ChIP-Seq ChIP-qPCR	Wiz	eLife	2016	Luke Isbel <i>et al.</i> Wiz binds active promoters and CTCF-binding sites and is required for normal behaviour in the mouse. <i>Elife.</i> 5: e15082
ChIP-Seq	BRD4, H3K27ac, H3K4me1, H3K27me3, HLX, LHX2, LMX1A	Nature	2016	Charles Y. Lin <i>et al.</i> Active medulloblastoma enhancers reveal subgroup specific cellular origins. <i>Nature.</i> 530: 57
ChIP-qPCR	H3K9me3 H3 pan-acetyl	Nucleic Acids Research	2016	Xiaoyu Chen <i>et al.</i> Probing the impact of chromatin conformation on genome editing tools. <i>Nucleic Acids Res.</i> 44: 6482
ChIP-Seq ChIP-qPCR	H3K9Ac	Stem Cells Translational Medicine	2016	Dalia Ali <i>et al.</i> Epigenetic library screen identifies abexinostat as novel regulator of adipocytic and osteoblastic differentiation of human skeletal (mesenchymal) stem cells. <i>Stem Cells Transl Med.</i> 5: 1036
ChIP-Seq	H3K27me3	Cancer Research	2016	John P. Thomson <i>et al.</i> Loss of Tet1 associated 5-hydroxymethylcytosine is concomitant with aberrant promoter hypermethylation in liver cancer. <i>Cancer Res.</i> Advance online publication doi: 10.1158/0008-5472.CAN-15-1910
ChIP-Seq, ChIP-qPCR	H3K36me3	Oncogene	2016	Thai H. Ho <i>et al.</i> High-resolution profiling of histone h3 lysine 36 trimethylation in metastatic renal cell carcinoma. <i>Oncogene.</i> 35: 1565
ChIP-Seq	HDAC1, HDAC2, GATA2	PLoS One	2016	Jeffrey R. Shearstone <i>et al.</i> Chemical Inhibition of histone deacetylases 1 and 2 Induces fetal hemoglobin through activation of GATA2. <i>PLoS One.</i> 11: e0153767
ChIP-Seq	LL-37 (antimicrobial peptide)	Journal of Cancer	2016	Mindy Munoz <i>et al.</i> Antimicrobial peptide LL-37 participates in the transcriptional regulation of melanoma cells. <i>J Cancer.</i> 26: 2341
ChIP-Seq ChIP-qPCR	p53	Molecular Oncology	2016	Cheryl Chan <i>et al.</i> Global re-wiring of p53 transcription regulation by the hepatitis B virus X protein. <i>Mol Oncol.</i> 10: 1183
ChIP-Seq	PLZF, SALL4	Development	2016	Dawn L. Lovelace <i>et al.</i> The regulatory repertoire of PLZF and SALL4 in undifferentiated spermatogonia. <i>Development.</i> Advance online publication doi: 10.1242/dev.132761
ChIP-Seq	PLZF	PLoS Genetics	2016	Ramakrishna Kommagani <i>et al.</i> The Promyelocytic leukemia zinc finger transcription factor is critical for human endometrial stromal cell decidualization. <i>PLoS Genet.</i> 12: e1005937
ChIP libraries, sequencing and analysis	H3K27me3	Developmental Dynamics	2016	Oyvind Dahle <i>et al.</i> Inhibiting smad2/3 signaling in pluripotent mouse embryonic stem cells enhances endoderm formation by increasing transcriptional priming of lineage-specifying target genes. <i>Dev Dyn.</i> 245: 807.
ChIP-Seq	NR4A1	PLoS One	2016	Ryan P. Duren <i>et al.</i> Genome wide mapping of NR4A binding reveals cooperativity with ETS factors to promote epigenetic activation of distal enhancers in acute myeloid leukemia cells. <i>PLoS One.</i> 11: e0150450

Technique	Target	Journal	Year	Reference
ChIP-Seq	MTA1	Oncotarget	2016	Swati Dhar <i>et al.</i> Dietary pterostilbene is a novel MTA1-targeted chemopreventive and therapeutic agent in prostate cancer. <i>Oncotarget</i> . Advance online publication doi: 10.18632/oncotarget.7841
ChIP-Seq	Myc-tagged Pet-1	The Journal of Neuroscience	2016	Steven C. Wyler <i>et al.</i> Pet-1 switches transcriptional targets postnatally to regulate maturation of serotonin neuron excitability. <i>J Neurosci</i> . 36:1758
ChIP-Seq	EGR2	Nature Communications	2015	Tomohisa Okamura <i>et al.</i> TGF- $\beta$ 3-expressing CD4 <sup>+</sup> CD25 <sup>-</sup> LAG3 <sup>+</sup> regulatory T cells control humoral immune responses. <i>Nat Commun</i> . 6: 6329
ChIP-Seq	HA-tagged Shox2	Development	2015	Wenduo Ye <i>et al.</i> A common Shox-2-Nkx2-5 antagonistic mechanism primes the pacemaker cell fate in the pulmonary vein myocardium and sinoatrial node. <i>Development</i> . 142: 2521
ChIP-Seq	BA180/PBRM1	Molecular Cell	2015	Bokai Zhu <i>et al.</i> Coactivator-dependent oscillation of chromatin accessibility dictates circadian gene amplitude via REV-ERB loading. <i>Mol Cell</i> . 60: 769
ChIP-qPCR	EZH2, SUZ12, H3K27me3, RNA Pol II	The EMBO Journal	2015	Stephen G. Dann <i>et al.</i> Reciprocal regulation of amino acid import and epigenetic state through Lat1 and EZH2. <i>EMBO J</i> . 34: 1773
ChIP-Seq	CCND1 (Cyclin D1)	Oncotarget	2015	Mathew C. Casimiro <i>et al.</i> Kinase-independent role of cyclin D1 in chromosomal instability and mammary tumorigenesis. <i>Oncotarget</i> . 6: 8524
ChIP-Seq	TRIM33	PLoS Genetics	2015	Luke Isbel <i>et al.</i> Trim33 binds and silences a class of young endogenous retroviruses in the mouse testis; a novel component of the arms race between retrotransposons and the host genome. <i>PLoS Genetics</i> . 11: e1005693
ChIP-Seq	Progesterone Receptor	Cell Reports	2015	Vincent J. Lynch <i>et al.</i> Ancient transposable elements transformed the uterine regulatory landscape and transcriptome during the evolution of mammalian pregnancy. <i>Cell Rep</i> . 10: 551
ChIP-Seq	EGR1	PLoS One	2015	Anthony A. Portale <i>et al.</i> Characterization of FGF23-dependent Egr-1 cistrome in the mouse renal proximal tubule. <i>PLoS One</i> . 10: e0142924
ChIP-Seq	FOXM1	Breast Cancer Research and Treatment	2015	Christina Yau <i>et al.</i> FOXM1 cistrome predicts breast cancer metastatic outcome better than FOXM1 expression levels or tumor proliferation index. <i>Breast Cancer Res Treat</i> . 154: 23
ChIP-Seq	BRD4	Molecular Cancer Therapy	2015	Ryan Lenhart <i>et al.</i> Sensitivity of small cell lung cancer to BET inhibition is mediated by regulation of ASCL1 gene expression. <i>Mol Cancer Ther</i> . 14: 2167
Next-Gen Bisulfite Seq	—	The Journal of Clinical Investigation	2015	Coralie Hoareau-Aveilla <i>et al.</i> Reversal of microRNA-150 silencing disadvantages crizotinib-resistant NPM-ALK(+) cell growth. <i>J Clin Invest</i> . 125:3505
ChIP-Seq ChIP-qPCR	b-Arrestin-1	Cancer Research	2015	Smitha Pillai <i>et al.</i> b-arrestin-1 mediates nicotine-induced metastasis through E2F1 target genes that modulate epithelial-mesenchymal transition. <i>Cancer Res</i> . 75: 1009
ChIP-Seq	H3K9ac, RNA Pol II, TFIIIB	Circulation: Heart Failure	2015	Danish Sayed <i>et al.</i> Acute targeting of general transcription factor IIB restricts cardiac hypertrophy via selective inhibition of gene transcription. <i>Circ Heart Fail</i> . 8: 138

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K4me3	Journal of the National Cancer Institute	2015	Cristian Taccioli <i>et al.</i> Repression of esophageal neoplasia and inflammatory signaling by anti-miR-31 delivery in vivo. <i>J Natl Cancer Inst.</i> 107: djv220
ChIP-Seq	FLAG-tagged Twist2	American Journal of Human Genetics	2015	Shannon Marchegiani <i>et al.</i> Recurrent mutations in the basic domain of TWIST2 cause ablepharon macrostomia and barber-say syndromes. <i>Am J Hum Genet.</i> 97: 99
ChIP-Seq	LHX6	Human Molecular Genetics	2015	Jeffry M. Cesario <i>et al.</i> Lhx6 and Lhx8 promote palate development through negative regulation of a cell cycle inhibitor gene, p57 <sup>Kip2</sup> . <i>Hum Mol Genet.</i> 24: 5024
ChIP-Seq	SOX9	Nucleic Acids Research	2015	Zhongcheng Shi <i>et al.</i> Context-specific role of SOX9 in NF-Y mediated gene regulation in colorectal cancer cells. <i>Nucleic Acids Res.</i> 43: 6257
ChIP-Seq	H3K27me3	Nucleic Acids Research	2015	Kyung Hyun Yoo <i>et al.</i> Loss of EZH2 results in precocious mammary gland development and activation of STAT5-dependent genes. <i>Nucleic Acids Res.</i> 43: 8774
ChIP-Seq	H3K4me1	BMC Biology	2015	Sara K. Harten <i>et al.</i> The recently identified modifier of murine metastable epialleles, Rearranged L-Myc Fusion, is involved in maintaining epigenetic marks at CpG island shores and enhancers. <i>BMC Biol.</i> 13: 21
ChIP-Seq	ZNF384	Molecular Endocrinology	2015	Paul Childress <i>et al.</i> Genome-wide mapping and interrogation of the Nmp4 anti-anabolic bone axis. <i>Mol Endocrinol.</i> 29: 1269
ChIP-Seq	FOXO1, RNA Pol II	Molecular Endocrinology	2015	Yasmin M. Vasquez <i>et al.</i> FOXO1 is Required for Binding of PR on IRF4, Novel Transcriptional Regulator of Endometrial Stromal Decidualization. <i>Mol Endocrinol.</i> 29: 421
ChIP-Seq	Progesterone Receptor, FOSL2	Endocrinology	2015	Erik C. Mazur <i>et al.</i> Progesterone receptor transcriptome and cistrome in decidualized human endometrial stromal cells. <i>Endocrinology.</i> 156: 2239
ChIP-Seq		Immunity	2015	Chrysothemis C. Brown <i>et al.</i> Retinoic acid is essential for Th1 cell lineage stability and prevents transition to a Th17 cell program. <i>Immunity.</i> 42: 1
ChIP-Seq	SIX1	Cancer Cell	2015	Jenny Wegert <i>et al.</i> Mutations in the SIX1/2 pathway and the DROSHA/DGCR8 miRNA microprocessor complex underlie high-risk blastemal type Wilms tumors. <i>Cancer Cell.</i> 27: 298
ChIP-Seq	H3K4me3, H3K9me3	Epigenetics	2015	Nioka C. Chisholm <i>et al.</i> Histone methylation patterns in astrocytes are influenced by age following ischemia. <i>Epigenetics.</i> 10: 142
ChIP-Seq Data Analysis	KDM3A	PLoS Biology	2015	Mo-bin Cheng <i>et al.</i> Specific phosphorylation of histone demethylase KDM3A determines target gene expression in response to heat shock. <i>PLOS Biol.</i> 12: e1002026
ChIP-Seq	SRC1	Molecular Endocrinology	2014	Mounia Tannour-Louet <i>et al.</i> Hepatic SRC-1 activity orchestrates transcriptional circuitries of amino acid pathways with potential relevance for human metabolic pathogenesis. <i>Mol Endocrinol.</i> 28: 1707
ChIP-Seq	Fish H3K4me3	Molecular Ecology Resources	2014	Claudius F. Kratochwil <i>et al.</i> Mapping active promoters by ChIP-seq profiling of H3K4me3 in cichlid fish – a first step to uncover cis-regulatory elements in ecological model teleosts. <i>Mol Ecol Resour.</i> Advance online publication. doi: 10.1111/1755-0998

Technique	Target	Journal	Year	Reference
ChIP-Seq	GATA2, FOXA1, p300, CBP, SRC1, SRC2, SRC3, Androgen Receptor	Proceedings of the National Academy of Sciences	2014	Bin He <i>et al.</i> GATA2 facilitates steroid receptor coactivator recruitment to the androgen receptor complex. <i>Proc Natl Acad Sci.</i> 111: 18261
Next-Gen Bisulfite Seq	—	Cell Reports	2014	Deepak Kumar <i>et al.</i> Fibroblast growth factor maintains chondrogenic potential of limb bud mesenchymal cells by modulating DNMT3A recruitment. <i>Cell Rep.</i> 8: 1419
ChIP-Seq	PPAR $\gamma$	Biology of Reproduction	2014	Kelsey E. Brooks <i>et al.</i> Peroxisome proliferator activator receptor gamma (PPARG) regulates conceptus elongation in sheep. <i>Biol Reprod.</i> 92: 42.
ChIP-Seq,	PXR, p300, H3K4me1, H3K27ac	PLoS Genetics	2014	Robin P. Smith <i>et al.</i> Genome-wide discovery of drug-dependent human liver regulatory elements. <i>PLOS Genet.</i> 10: e1004648
ChIP-Seq	H3K9me3, H3K36me3, H3K27me3	Genome Research	2014	Goran Kungulovski <i>et al.</i> Application of histone modification-specific interaction domains as an alternative to antibodies. <i>Genome Res.</i> 24: 1842
ChIP-Seq ChIP-qPCR	SRF	Physiological Genomics	2014	Sharolyn V. Kawakami-Schulz. Serum response factor: positive and negative regulation of an epithelial gene expression network in the destrin mutant cornea. <i>Physiol Genomics.</i> 46: 277
ChIP-Seq ChIP-qPCR	SRC3	The Journal of Clinical Investigation	2014	Jun Qin <i>et al.</i> Androgen deprivation–induced NCoA2 promotes metastatic and castration-resistant prostate cancer. <i>J Clin Invest.</i> 124: 5013
ChIP-Seq	H3K27ac	Journal of Biological Chemistry	2014	Andre Landin Malt <i>et al.</i> Identification of a face enhancer reveals direct regulation of LIM homeobox 8 (Lhx8) by wingless-int (WNT)/b-catenin Signaling. <i>J Biol Chem.</i> 289: 30289
ChIP-Seq ChIP-qPCR	ZBTB20	Journal of Medical Genetics	2014	Malene B. Rasmussen <i>et al.</i> Neurodevelopmental disorders associated with dosage imbalance of ZBTB20 correlate with the morbidity spectrum of ZBTB20 candidate target genes. <i>J Med Genet.</i> 51: 605
ChIP-Seq	H3K9Ac, H3K27ac	Nature	2014	Paul A. Northcott <i>et al.</i> Enhancer hijacking activates GF1 family oncogenes in medulloblastoma. <i>Nature.</i> 511: 428
ChIP-Seq	H3K4me3, H3K27me3	Nature	2014	Volker Hovestadt <i>et al.</i> Decoding the regulatory landscape of medulloblastoma using DNA methylation sequencing. <i>Nature.</i> 510: 537
ChIP-qPCR	EZH2, Suz12, H3K27me3	Molecular Cancer Therapy	2014	Sarah K. Knutson <i>et al.</i> Selective Inhibition of EZH2 by EPZ-6438 Leads to Potent Antitumor Activity in EZH2 Mutant Non-Hodgkin Lymphoma. <i>Mol Cancer Ther.</i> 13: 842
ChIP-Seq	RAR-a, RAR-b, RXR-a	American Journal of Physiology – Gastrointestinal and Liver Physiology	2014	Yuqi He <i>et al.</i> Biological functional annotation of retinoic acid alpha and beta in mouse liver based on genome-wide binding. <i>Am J Physiol Gastrointest Liver Physiol.</i> 307: G205
ChIP-Seq	H3K4me3, H3K27me3	Nature Communications	2014	Brian C. Belyea <i>et al.</i> Identification of renin progenitors in the mouse bone marrow that give rise to B-cell leukaemia. <i>Nat Commun.</i> 5: 3273
ChIP-Seq	Estrogen Receptor	Molecular Endocrinology	2014	Sylvia C. Hewitt <i>et al.</i> Novel DNA motif binding activity observed In vivo with an estrogen receptor a mutant mouse. <i>Mol Endocrinol.</i> 26: 899

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27me3	Bioinformatics	2014	Yanxiao Zhang <i>et al.</i> PePr: A peak-calling prioritization pipeline to identify consistent or differential peaks from replicated ChIP-Seq data. <i>Bioinformatics</i> . 30: 2568
ChIP-Seq	H3K36me3	Genome Research	2014	Jeremy M. Simon <i>et al.</i> Variation in chromatin accessibility in human kidney cancer links H3K36 methyltransferase loss with widespread RNA processing defects. <i>Genome Res.</i> 24: 241
ChIP-Seq	SRC-2	Cell Reports	2014	Erin Stashi <i>et al.</i> SRC-2 Is an Essential Coactivator for Orchestrating Metabolism and Circadian Rhythm. <i>Cell Rep.</i> 6: 663
Next-Gen Bisulfite Seq	—	Nature Neuroscience	2014	Brian G. Dias <i>et al.</i> Parental olfactory experience influences behavior and neural structure in subsequent generations. <i>Nat Neurosci.</i> 17: 89
ChIP-Seq ChIP-qPCR	Zbtb20	Cerebral Cortex	2014	Jakob V. Nielsen <i>et al.</i> Zbtb20 Defines a Hippocampal Neuronal Identity Through Direct Repression of Genes That Control Projection Neuron Development in the Isocortex. <i>Cereb Cortex.</i> 24: 1216.
ChIP-qPCR	H3K9me3, H3	Nature Immunology	2013	Patrick M. Gubser <i>et al.</i> Rapid effector function of memory CD8+ T cells requires an immediate-early glycolytic switch. <i>Nat Immunol.</i> 14: 1064
ChIP-Seq	RNA Pol II	PLoS One	2013	Jonathan P. Riley <i>et al.</i> PARP-14 Binds Specific DNA Sequences to Promote Th2 Cell Gene Expression. <i>PLoS One.</i> 8: e83127
Next-Gen Bisulfite Seq	—	Clinical Cancer Research	2013	David S. Shames <i>et al.</i> Loss of NAPRT1 Expression by Tumor-Specific Promoter Methylation Provides a Novel Predictive Biomarker for NAMPT Inhibitors. <i>Clin Cancer Res.</i> 19: 6912
ChIP-Seq	Cyclin D1	Cancer Research	2013	Xiaoming Ju <i>et al.</i> Identification of a cyclin D1 network in prostate cancer that antagonizes epithelial-mesenchymal restraint. <i>Cancer Res.</i> 74: 508
ChIP-Seq	H3K27me3	Cancer Cell	2013	Sebastian Bender <i>et al.</i> Reduced H3K27me3 and DNA Hypomethylation Are Major Drivers of Gene Expression in K27M Mutant Pediatric High-Grade Gliomas. <i>Cancer Cell.</i> 24: 660
ChIP-Seq	ASXL1	Journal of Experimental Medicine	2013	Omar Abdel-Wahab <i>et al.</i> Deletion of Asxl1 results in myelodysplasia and severe developmental defects in vivo. <i>J Exp Med.</i> 210: 2641
ChIP-Seq	COUP-TFII (NR2F2)	Molecular Endocrinology	2013	Xilong Li <i>et al.</i> COUP-TFII Regulates Human Endometrial Stromal Genes Involved in Inflammation. <i>Mol Endocrinol.</i> 27: 2041
ChIP-Seq	Gfi1	Nature Immunology	2013	Chauncey J. Spooner <i>et al.</i> Specification of type 2 innate lymphocytes by the transcriptional determinant Gfi1. <i>Nat Immunol.</i> 14: 1229
ChIP-Seq	RNA Pol II	Journal of Allergy and Clinical Immunology	2013	Purvi Mehrotra <i>et al.</i> PARP-14 and its enzyme activity regulates Th2 differentiation and allergic airway disease. <i>J Allergy Clin Immunol.</i> 131: 52
ChIP-Seq	Progesterone Receptor	Molecular Endocrinology	2013	Ashlee R. Lain <i>et al.</i> Research Resource: Progesterone Receptor Targetome Underlying Mammary Gland Branching Morphogenesis. <i>Mol Endocrinol.</i> 27: 1743
ChIP-Seq	PPARα	Chemico-Biological Interactions	2013	Patrick D. McMullen <i>et al.</i> A map of the PPARα transcription regulatory network for primary human hepatocytes. <i>Chem Biol Interact.</i> 209: 104

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K36me3	Cancer Discovery	2013	Lynn Bjerke <i>et al.</i> Histone H3.3 Mutations Drive Pediatric Glioblastoma through Upregulation of MYCN. <i>Cancer Discov.</i> 3: 512
ChIP-Seq	p53	Physiological Genomics	2013	Yuwen Li <i>et al.</i> Genome-wide analysis of the p53 gene regulatory network in the developing mouse kidney. <i>Physiol Genomics.</i> 45: 948
ChIP-Seq	FOXA2	The FASEB Journal	2013	Justyna Filant <i>et al.</i> Integrated chromatin immunoprecipitation sequencing and microarray analysis identifies FOXA2 target genes in the glands of the mouse uterus. <i>FASEB J.</i> 28: 230
ChIP-Seq	p300	PLoS Genetics	2013	Aaron M. Wenger <i>et al.</i> The Enhancer Landscape during Early Neocortical Development Reveals Patterns of Dense Regulation and Co-option. <i>PLoS Genet.</i> 9: e1003728
ChIP-Seq	MYRF	PLoS Biology	2013	Helena Bujalka <i>et al.</i> MYRF is a membrane-associated transcription factor that autoproteolytically cleaves to directly activate myelin genes. <i>PLoS Biol.</i> 11: e1001625
ChIP-Seq ChIP-qPCR	RXR- $\alpha$ , RNA Pol II	PLoS One	2013	Astrid Kosters <i>et al.</i> Sexually Dimorphic Genome-Wide Binding of Retinoid X Receptor $\alpha$ (RXR $\alpha$ ) Determines Male-Female Differences in the Expression of Hepatic Lipid Processing Genes in Mice. <i>PLoS One.</i> 8: e71538
ChIP-qPCR	NF $\kappa$ B (p52 & p65)	Journal of Biological Chemistry	2013	Sarah L. Doyle <i>et al.</i> Nuclear factor $\kappa$ B2 p52 protein has a role in antiviral immunity through I $\kappa$ B kinase epsilon-dependent induction of Sp1 protein and interleukin 15. <i>J Biol Chem.</i> 288: 25066
ChIP-chip ChIP-qPCR	C/EBP $\beta$ , RNA Pol II	Journal of Biological Chemistry	2013	Hana Vakili <i>et al.</i> CCAAT-enhancer-binding protein b (C/EBP $\beta$ ) and downstream human placental growth hormone genes are targets for dysregulation in pregnancies complicated by maternal obesity. <i>J Biol Chem.</i> 288: 22849
ChIP-Seq	KDM2B	The Journal of Clinical Investigation	2013	Alexandros Tzatsos <i>et al.</i> KDM2B promotes pancreatic cancer via Polycomb-dependent and -independent transcriptional programs. <i>J Clin Invest.</i> 123: 727
ChIP-Seq	Vitamin D Receptor	BMC Medicine	2013	Adam E. Handel <i>et al.</i> Vitamin D receptor ChIP-seq in primary CD4 <sup>+</sup> cells: relationship to serum 25-hydroxyvitamin D levels and autoimmune disease. <i>BMC Med.</i> 11: 163
ChIP-Seq	COUP-TFII (NR2F2)	Developmental Cell	2013	San-pin Wu <i>et al.</i> Atrial Identity Is Determined by a COUP-TFII Regulatory Network. <i>Dev Cell.</i> 25: 417
ChIP-Seq	H3K4me3	PLoS One	2013	Alex Gutteridge <i>et al.</i> Novel Pancreatic Endocrine Maturation Pathways Identified by Genomic Profiling and Causal Reasoning. <i>PLoS One.</i> 8: e56024.
ChIP-Seq	Glucocorticoid Receptor	PLoS One	2012	Marie-José C. van Lierop <i>et al.</i> Org 214007-0: A Novel Non-Steroidal Selective Glucocorticoid Receptor Modulator with Full Anti- Inflammatory Properties and Improved Therapeutic Index. <i>PLoS One.</i> 7: e48385
ChIP-Seq	KDM1A (LSD1), H3K4me2	Molecular and Cellular Biology	2012	Venugopalan D. Nair <i>et al.</i> Involvement of Histone Demethylase LSD1 in Short-Time-Scale Gene Expression Changes during Cell Cycle Progression in Embryonic Stem Cells. <i>Mol Cell Biol.</i> 32: 4861
ChIP-Seq	EZH2	Nature	2012	Michael T. McCabe <i>et al.</i> EZH2 inhibition as a therapeutic strategy for lymphoma with EZH2-activating mutations. <i>Nature.</i> 492: 108



Technique	Target	Journal	Year	Reference
Data Analysis	pTyr37 H2B	Nature Structural and Molecular Biology	2012	Kiran Mahajan <i>et al.</i> H2B Tyr37 phosphorylation suppresses expression of replication-dependent core histone genes. <i>Nat Struct Mol Biol.</i> 19: 930
ChIP-chip	MYBL2 (B-MYB)	PLoS One	2012	Ming Zhan <i>et al.</i> The B-MYB Transcriptional Network Guides Cell Cycle Progression and Fate Decisions to Sustain Self-Renewal and the Identity of Pluripotent Stem Cells. <i>PLoS One.</i> 7: e42350
ChIP-Seq	FLAG-BAP1, OGT, HCF1	Science	2012	Anwasha Dey <i>et al.</i> Loss of the Tumor Suppressor BAP1 Causes Myeloid Transformation. <i>Science.</i> 337: 1541
ChIP-Seq	RNA Pol II	Journal of Biological Chemistry	2012	Danish Sayed <i>et al.</i> Transcriptional regulation patterns revealed by high-resolution chromatin immunoprecipitation during cardiac hypertrophy. <i>J Biol Chem.</i> 288: 2546
ChIP-Seq	Androgen Receptor	BMC Genomics	2012	Zhou Zhu <i>et al.</i> Dose-dependent effects of small-molecule antagonists on the genomic landscape of androgen receptor binding. <i>BMC Genomics.</i> 13: 355
ChIP-Seq	Progesterone Receptor	Molecular Endocrinology	2012	Cory A. Rubel <i>et al.</i> Genome-Wide Profiling of Progesterone Receptor Binding in the Mouse Uterus. <i>Mol Endocrinol.</i> 26: 1428
ChIP-Seq	Estrogen Receptor, RNA Pol II	Molecular Endocrinology	2012	Sylvia C. Hewitt <i>et al.</i> Whole-Genome Estrogen Receptor a Binding in Mouse Uterine Tissue Revealed by ChIP-Seq. <i>Mol Endocrinol.</i> 26: 887
ChIP-qPCR	PDX1, TCF3	Chemistry & Biology	2012	Alice Kiselyuk <i>et al.</i> HNF4a antagonists discovered by a high-throughput screen for modulators of the human insulin promoter. <i>Chem Biol.</i> 19: 806
ChIP-chip	RORa	PLoS One	2012	Yongjun Wang <i>et al.</i> Regulation of p53 Stability and Apoptosis by a ROR Agonist. <i>PLoS One.</i> 7: e34921
ChIP-Seq	Nkx3.1	The Journal of Clinical Investigation	2012	Philip D. Anderson <i>et al.</i> Nkx3.1 and Myc crossregulate shared target genes in mouse and human prostate tumorigenesis. <i>J Clin Invest.</i> 122: 1907
ChIP-Seq	RNA Pol II	Cardiovascular Research	2012	Mingyue Han <i>et al.</i> GATA4 expression is primarily regulated via a miR-26b-dependent post-transcriptional mechanism during cardiac hypertrophy. <i>Cardiovasc Res.</i> 93: 645
ChIP-Seq	FXR (Farnesoid X Receptor)	The American Journal of Physiology: Gastrointestinal and Liver Physiology	2012	Julia Yue Cui <i>et al.</i> Bile acids via FXR initiate the expression of major transporters involved in the enterohepatic circulation of bile acids in newborn mice. <i>Am J Physiol Gastrointest Liver Physiol.</i> 302: G979
ChIP-Seq	Progesterone Receptor	PLoS One	2012	Ping Yin <i>et al.</i> Genome-wide progesterone receptor binding: cell type-specific and shared mechanisms in T47D breast cancer cells and primary leiomyoma cells. <i>PLoS One.</i> 7: e29021
ChIP-Seq	FLAG-CCND1 (Cyclin D1)	The Journal of Clinical Investigation	2012	Mathew C. Casimiro <i>et al.</i> ChIP sequencing of cyclin D1 reveals a transcriptional role in chromosomal instability in mice. <i>J Clin Invest.</i> 122: 833
ChIP-Seq	RBPJ	Stem Cells	2012	Yaochen Li <i>et al.</i> Genome-wide Analysis of N1ICD/RBPJ Targets In Vivo Reveals Direct Transcriptional Regulation of Wnt, SHH, and Hippo Pathway Effectors by Notch1. <i>Stem Cells.</i> 30: 741

Technique	Target	Journal	Year	Reference
ChIP-chip	Progesterone Receptor	The FASEB Journal	2011	Heather L. Franco <i>et al.</i> Epithelial progesterone receptor exhibits pleiotropic roles in uterine development and function. <i>FASEB J.</i> 26: 1218
ChIP-Seq	Msgn1	Nature Communication	2011	Ravindra B. Chalamalasetty <i>et al.</i> The Wnt3a/b-catenin target gene Mesogenin1 controls the segmentation clock by activating a Notch signalling program. <i>Nat Commun.</i> 2: 390
ChIP-qPCR	IRF4, Myc	British Journal of Haematology	2011	Antonia Lopez-Girona <i>et al.</i> Lenalidomide downregulates the cell survival factor, interferon regulatory factor-4, providing a potential mechanistic link for predicting response. <i>Br J Haematol.</i> 154: 325
ChIP-chip	IRF4	Allergy	2011	Barrenas S. Bruhn <i>et al.</i> Increased expression of IRF4 and ETS1 in CD4+ cells from patients with intermittent allergic rhinitis. <i>Allergy.</i> 67: 33
ChIP-Seq	COBRA1 (Nelf-b)	Journal of Biological Chemistry	2011	Jianlong Sun <i>et al.</i> Genetic and Genomic Analyses of RNA Polymerase II-pausing Factor in Regulation of Mammalian Transcription and Cell Growth. <i>J Biol Chem.</i> 286: 36248
ChIP-Seq	Foxp3	Nucleic Acids Research	2011	Fabian Birzele <i>et al.</i> Next-generation insights into regulatory T cells: expression profiling and FoxP3 occupancy in Human. <i>Nucleic Acids Res.</i> 39: 7946
ChIP-qPCR	BRD4	Proceedings of the National Academy of Sciences	2011	Jennifer A. Mertz <i>et al.</i> Targeting MYC dependence in cancer by inhibiting BET bromodomains. <i>Proc Natl Acad Sci.</i> 108: 16669
ChIP-chip	C/EBPb	Journal of Molecular Endocrinology	2011	Aristides Lytras <i>et al.</i> Identification of functional CCAAT/ enhancer-binding protein and Ets protein binding sites in the human chorionic somatomammotropin enhancer sequences. <i>J Mol Endocrinol.</i> 47: 179
ChIP-qPCR	Progesterone Receptor	Biochemical Pharmacology	2011	Matthew R. Yudit <i>et al.</i> Discovery of a novel mechanism of steroid receptor antagonism: WAY-255348 modulates progesterone receptor cellular localization and promoter interactions. <i>Biochem Pharmacol.</i> 82: 1709
ChIP-chip	RNA Pol II, H3K9Ac	Journal of Biological Chemistry	2011	Hong Hao <i>et al.</i> The Transcription Factor Neural Retina Leucine Zipper (NRL) Controls Photoreceptor-specific Expression of Myocyte Enhancer Factor Mef2c from an Alternative Promoter. <i>J Biol Chem.</i> 286: 34893
ChIP-chip	DAF-12	PLoS Genetics	2011	Daniel Hochbaum <i>et al.</i> DAF-12 Regulates a Connected Network of Genes to Ensure Robust Developmental Decisions. <i>PLoS Genet.</i> 7: e1002179
ChIP-qPCR	CTCF	DNA and Cell Biology	2011	Yan Jin <i>et al.</i> Enhancer-Blocking Activity Is Associated with Hypersensitive Site V Sequences in the Human Growth Hormone Locus Control Region. <i>DNA Cell Biol.</i> 30: 995
Bisulfite-Seq	Bisulfite Sequencing	PLoS One	2011	Michela Deleidi <i>et al.</i> Oct4-Induced Reprogramming Is Required for Adult Brain Neural Stem Cell Differentiation into Midbrain Dopaminergic Neurons. <i>PLoS ONE.</i> 6: e19926
MeDIP-chip	5-methylcytosine (5-mC)	BMC Biology	2011	Genevieve Lavoie <i>et al.</i> PKC isoforms interact with and phosphorylate DNMT1. <i>BMC Biol.</i> 9: 31
ChIP-qPCR	AHR (Aryl Hydrocarbon Receptor)	Toxicology and Applied Pharmacology	2011	K. Nadira De Abrew <i>et al.</i> Regulation of Bach2 by the aryl hydrocarbon receptor as a mechanism for suppression of B-cell differentiation by 2,3,7,8-tetrachlorodibenzo-p-dioxin. <i>Toxicol Appl Pharmacol.</i> 252: 150

Technique	Target	Journal	Year	Reference
ChIP-qPCR	PPAR $\gamma$	Investigative Ophthalmology and Visual Science	2011	Gerard A. Rodrigues <i>et al.</i> Differential Effects of PPAR $\gamma$ Ligands on Oxidative Stress-Induced Death of Retinal Pigmented Epithelial Cells. <i>Invest Ophthalmol Vis Sci.</i> 52: 890
ChIP-chip	p73	Proceedings of the National Academy of Sciences	2011	Jennifer M. Rosenbluth <i>et al.</i> Differential regulation of the p73 cistrome by mammalian target of rapamycin reveals transcriptional programs of mesenchymal differentiation and tumorigenesis. <i>Proc Natl Acad Sci.</i> 108: 2076
ChIP-chip	LXR, RXR	Journal of Biological Chemistry	2011	Qi Shen <i>et al.</i> Liver X Receptor-Retinoid X Receptor (LXR-RXR) Heterodimer Cistrome Reveals Coordination of LXR and AP1 Signaling in Keratinocytes. <i>J Biol Chem.</i> 286: 14554
ChIP-qPCR	AHR (Aryl Hydrocarbon Receptor), MAF	Nature Immunology	2010	Lionel Apetoh <i>et al.</i> The aryl hydrocarbon receptor interacts with c-Maf to promote the differentiation of type 1 regulatory T cells induced by IL-27. <i>Nat Immunol.</i> 11: 854
ChIP-qPCR	p73	Molecular Cancer	2010	Kathryn G. Eby <i>et al.</i> ISG20L1 is a p53 family target gene that modulates genotoxic stress-induced autophagy. <i>Mol Cancer.</i> 9: 95
ChIP-Seq	Vitamin D Receptor	Genome Research	2010	Sreeram V. Ramagopalan <i>et al.</i> A ChIP-seq defined genome-wide map of vitamin D receptor binding: associations with disease and evolution. <i>Genome Res.</i> 20: 1352
ChIP-chip	O-GlcNAc, RNA Pol II	Proceedings of the National Academy of Sciences	2010	Dona C. Love <i>et al.</i> Dynamic O-GlcNAc cycling at promoters of <i>Caenorhabditis elegans</i> genes regulating longevity, stress, and immunity. <i>Proc Natl Acad Sci.</i> 107: 7413
ChIP-Seq	PXR (Pregnane X Receptor)	Toxicological Sciences	2010	Julia Yue Cui <i>et al.</i> Genetic and Epigenetic Regulation and Expression Signatures of Glutathione S-Transferases in Developing Mouse Liver. <i>Toxicol Sci.</i> 116: 32
ChIP-Seq	SRC3, RNA Pol II	Molecular Endocrinology	2010	Rainer B. Lanz <i>et al.</i> Global Characterization of Transcriptional Impact of the SRC-3 Coregulator. <i>Mol Endocrinol.</i> 24: 859
ChIP-qPCR	Maf1, Rpc39, mTOR, Brf1, Raptor	Journal of Biological Chemistry	2010	Boris Shor <i>et al.</i> Requirement of the mTOR Kinase for the Regulation of Maf1 Phosphorylation and Control of RNA Polymerase III-dependent Transcription in Cancer Cells. <i>J Biol Chem.</i> 285: 15380
ChIP-chip	AHR (Aryl Hydrocarbon Receptor)	Toxicological Sciences	2010	K. Nadira De Abrew <i>et al.</i> An Integrated Genomic Analysis of Aryl Hydrocarbon Receptor-Mediated Inhibition of B-Cell Differentiation. <i>Toxicol Sci.</i> 118: 454
ChIP-qPCR	Myc-tagged-Pet-1	Nature Neuroscience	2010	Chen Liu <i>et al.</i> Pet-1 is required across different stages of life to regulate serotonergic function. <i>Nat Neurosci.</i> 13: 1190
ChIP-chip	Androgen Receptor	Molecular Endocrinology	2010	Anastasia Wyce <i>et al.</i> The Androgen Receptor Modulates Expression of Genes with Critical Roles in Muscle Development and Function. <i>Mol Endocrinol.</i> 24: 1665
ChIP-qPCR	NF $\kappa$ B (p50 and p65)	Cardiovascular Research	2010	Alina G. Sofronescu <i>et al.</i> FGF-16 is a target for adrenergic stimulation through NF- $\kappa$ B activation in postnatal cardiac cells and adult mouse heart. <i>Cardiovasc Res.</i> 87: 102
ChIP-Seq	PXR (Pregnane X Receptor)	Nucleic Acids Research	2010	Julia Yue Cui <i>et al.</i> ChIPing the cistrome of PXR in mouse liver. <i>Nucleic Acids Res.</i> 38: 7943
ChIP-chip	RNA Pol II	Molecular Vision	2010	Padmaja Tummala <i>et al.</i> Temporal ChIP-on-Chip of RNA-Polymerase-II to detect novel gene activation events during photoreceptor maturation. <i>Mol Vis.</i> 16: 252

Technique	Target	Journal	Year	Reference
ChIP-chip	STAT4	The Journal of Immunology	2009	Seth R. Good <i>et al.</i> Temporal Induction Pattern of STAT4 Target Genes Defines Potential for Th1 Lineage-Specific Programming. <i>J Immunol.</i> 183: 3839
ChIP-Seq	p73	Molecular and Cellular Biology	2008	Jennifer M. Rosenbluth <i>et al.</i> A Gene Signature-Based Approach Identifies mTOR as a Regulator of p73. <i>Mol Cell Biol.</i> 28: 5951
ChIP-qPCR	AHR (Aryl Hydrocarbon Receptor)	Nature	2008	Francisco J. Quintana <i>et al.</i> Control of Treg and TH17 cell differentiation by the aryl hydrocarbon receptor. <i>Nature.</i> 453: 65