

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.

Technique	Target	Journal	Year	Reference
ATAC-Seq	ATAC-seq	Nature	2020	Biraj Mahato et al. 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 et al. 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 et al. 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 et al. MRG15 orchestrates rhythmic epigenomic remodelling and controls hepatic lipid metabolism. <i>Nat Metab</i> . 2020 doi: 10.1038/s42255-020-0203-z
HiC	NA	Journal of Biological Chemistry	2020	Sylvia Hewitt et al. Estrogen receptor α (ER α)-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	NA	PNAS	2020	Kruttika Bhat et al. 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
ATAC-Seq, ChIP-Seq	N-MYC, C-MYC, H3K4me1, H3K4me3, H3K27ac, and H3K27me3	Scientific Data	2020	Kristen Upton et al. Epigenomic profiling of neuroblastoma cell lines. <i>Sci Data</i> . 2020 doi: 10.6084/m9.figshare.11993883
ChIP-Seq	H3K27ac and AFF4	Communications Biology	2020	Yujing Gao et al. 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, and RNA Pol II	Experimental Cell Research	2020	Paul-Joseph Aspuria et al. 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 et al. 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	NA	Cancer Discovery	2020	Di Zhao et al. 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	NA	Diabetes	2020	Marion de Toledo et al. Lamin C counteracts glucose intolerance in aging, obesity and diabetes. <i>Diabetes</i> . 2020 doi: 10.2337/db19-0377.
ChIP-Seq	N-MYC and H3K27ac	Neoplasia	2020	Dongdong Chen et al. 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 et al. CAMKV Is a Candidate Immunotherapeutic Target in MYCN Amplified Neuroblastoma. <i>Front. Oncol</i> . 2020 doi: 10.3389/fonc.2020.00302

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K4me3, H3K9me3, H3K27me3, H3R8me2a, H4K20me3	Theranostics	2020	Qian Zhang et al. 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 et al. 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 et al. 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	NA	Cell Reports	2019	Banushree Kumar et al. 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	NA	The EMBO Journal	2019	Kathryn A. Jacobs et al. 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 et al. Therapeutic ligands antagonize estrogen receptor function by impairing its mobility. <i>Cell</i> . 178: 15
ATAC-Seq, ChIP-Seq	H3K4me3, H3K27Ac, H3K27me3	Nature	2019	Nina Oberbeck et al. The RIPK4–IRF6 signalling axis safeguards epidermal differentiation and barrier function. <i>Nature</i> . 574: 249
ATAC-Seq	NA	Cancer Cell	2019	Ana P. Gomes et al. 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 et al. 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 et al. 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 et al. 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 et al. 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 et al. 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 et al. 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 et al. Nucleome dynamics during retinal development. <i>Neuron</i> . 104:1
ChIP-qPCR	NRF2	Nature Immunology	2019	Anne-Valerie Burgener et al. 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 et al. 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"

Technique	Target	Journal	Year	Reference
ChIP-Seq	H4 panAc	Cell Metabolism	2019	Ramon C. Sun et al. 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 et al. 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 et al. The COMPASS Family Protein ASH2L Mediates Corticogenesis via Transcriptional Regulation of Wnt Signaling. <i>Cell Reports.</i> 28: 698
Mod Spec	NA	Nature Genetics	2019	Sara Martire et al. 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 et al. Long-read sequencing unveils IGH-DUX4 translocation into the silenced IGH allele in B-cell acute lymphoblastic leukemia. <i>Nat. Commun.</i> 10: 2789
ChIP-Seq	SALL3, DNMT3B	Nature Communications	2019	Takuya Kuroda et al. 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 et al. 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 et al. 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	NA	Haematologica	2019	James Ropa et al. 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 et al. 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 et al. 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 et al. 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 et al. 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 et al. 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 et al. Tissue-specific progesterone receptor-chromatin binding and the regulation of progesterone-dependent gene expression. <i>Sci Rep.</i> 19: 9
Next-Gen Bisulfite-Seq	NA	Plant Biotechnology Journal	2019	Aurine Verkest et al. 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

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K9me3	Journal of Experimental Medicine	2019	Jun Hyung Sin et al. The epigenetic regulator ATF7ip inhibits IL2 expression, regulating Th17 responses. <i>J. Exp Med.</i> 216: 2024
ChIP-Seq	H3K27Ac	Cancer Research	2019	Joeseph Castillo et al. 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 et al. Induction of a cell-restricted Gc in dedifferentiating β cells contributes to stress-induced β -cell dysfunction. <i>J Clin Invest.</i> 5: 128351
ChIP-Seq	SMC1a, H3K4me1, H3K4me3, H3K27Ac	Journal of Biological Chemistry	2019	Sylvia C. Hewitt et al. 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
ChIP-Seq, ChIP-qPCR	CBP, p300, H3K27Ac	Cell Reports	2019	Denise de Almeida Nagata et al. 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	Glucocorticoid Receptor	Journal of the American Heart Association	2019	Elena Severinova et al. Glucocorticoid Receptor-Binding and Transcriptome Signature in Cardiomyocytes. <i>J Am Heart Assoc.</i> 8: e011484.
ChIP-Seq	H4K20me1	Journal of Autoimmunity	2019	Keqi Fan et al. CRL4DCAF2is required for mature T-cell expansion via Aurora B-regulated proteasome activity. <i>J of Autoimmun.</i> 96:74
MeDIP-Seq	5-Methylcytosine	Epigenomes	2019	Liliana Ferreira et al. 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 et al. Interaction between hormone-sensitive lipase and ChREBP in fat cells controls insulin sensitivity. <i>Nat. Metab.</i> 1: 133
ChIP-Seq	H3K4me1, H3K-4me3, H3K27Ac	Cell Metabolism	2019	Liming Du et al. IGF-2 Preprograms Maturing Macrophages to Acquire Oxidative Phosphorylation-Dependent Antiinflammatory Properties. <i>Cell Metab.</i> 29: 1
ChIP-Seq	RNA Pol II, H3K27Ac	Journal of Neuro-Oncology	2019	Isabel Tegeder et al. Functional relevance of genes predicted to be affected by epigenetic alterations in atypical teratoid/rhabdoid tumors. <i>J. Neurooncol.</i> 141: 43
ChIP-Seq	EGR1, H3K27Ac	Journal of Steroid Biochemistry & Molecular Biology	2019	Maria Szwarc et al. Early growth response 1 transcriptionally primes the human endometrial stromal cell for decidualization. <i>J Steroid Biochem.</i> 189: 283
ChIP-Seq	H3K27me3, H3K-4me3	Acta Neuropathologica	2019	Andre Silveira et al. H3.3 K27M depletion increases differentiation and extends latency of diffuse intrinsic pontine glioma growth in vivo. <i>Acta Neuropathol.</i> 137: 1021
ChIP-Seq	RNA Pol II	Agronomy	2019	Sonja Klemme et al. Selection of Salicylic Acid Tolerant Epilines in Brassica napus. <i>Agronomy.</i> 9: 92
ChIP-Seq	LSD1, H3K4me2, H3K27Ac	Science Signaling	2019	Arnaud Augert et al. Targeting NOTCH activation in small cell lung cancer through LSD1 inhibition. <i>Sci.Signal.</i> 12: eaau2922
ChIP-qPCR	SRE, CEBP	Nature Communications	2019	S. Sarcar et al. Next-generation muscle-directed gene therapy by in silico vector design. <i>Nat Commun.</i> 10: 492
ChIP-Seq	H3K27Ac, H3K4me3, CTCF, NR4A3	Nature Communications	2019	Florian Haller et al. Enhancer hijacking activates oncogenic transcription factor NR4A3 in acinic cell carcinomas of the salivary glands. <i>Nat Commun.</i> 10: 336

Technique	Target	Journal	Year	Reference
ChIP-Seq	GFI1, LSD1	Nature Communications	2019	Catherine Lee et al. Lsd1 as a therapeutic target in Gfi1-activated medulloblastoma. <i>Nat Commun.</i> 10: 332
ChIP-Seq	H3K27me3, H3K4me1, H3K4me3, H3K9me3, H3K36me3, EZH2, SMARCB1, SUZ12, SMARCA4, REST	Cancer Cell	2019	Serap Erkek et al. 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 et al. Chromatin remodeling ATPase BRG1 and PTEN are synthetic lethal in prostate cancer. <i>J Clin Invest.</i> 129: 759
ChIP-qPCR	PPARg	International Journal of Obesity	2019	Laura Butruille et al. 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 (London)</i> doi: 10.1038/s41366-018-0310-z
ChIP-Seq	HA-tagged JP2NT, TBP, MEF2C	Science	2018	Ang Guo et al. 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 et al. A proteasome-resistant fragment of NIK mediates oncogenic NF-κB signaling in schwannomas. <i>Hum Mol Gen.</i> 28:572
ChIP-Seq	NRF2	Journal of Biological Chemistry	2018	Junsheng Fu et al. 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 et al. 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 et al. 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 Wang et al. 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	NA	Cell Reports	2018	Xiaolong Zhang et al. OX40 Costimulation inhibits Foxp3 expression and Treg induction via BATF3-dependent and independent mechanisms. <i>Cell Reports.</i> 24: 607
Mod Spec	NA	Acta Neuropathologica	2018	Leah Katz et al. Loss of histone H3K27me3 identifies a subset of meningiomas with increased risk of recurrence <i>Acta Neuropathol.</i> 135:955
RIME	TRPS1	Cell Reports	2018	Robert M. Witwicki et al. TRPS1 Is a lineage-specific transcriptional dependency in breast cancer. <i>Cell Reports.</i> 25: 1255
ChIP-Seq	H3K27Ac	Cell	2018	Xun Huang et al. 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 et al. Identification of therapeutic targets in rhabdomyosarcoma through integrated genomic, epigenomic, and proteomic analyses. <i>Cancer Cell.</i> 34: 1

Technique	Target	Journal	Year	Reference
ChIP-Seq	SOX17	Nature Communications	2018	Xiaoqiu Wang et al. SOX17 regulates uterine epithelial-stromal cross-talk acting via a distal enhancer upstream of <i>Ihh</i> . <i>Nat Commun.</i> 9: 4421
ChIP-Seq, RNA-Seq	NRF2, RNA Pol II	Nature Communications	2018	David Olagnier et al. 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 et al. 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 et al. Super-Enhancers maintain renin-expressing cell identity and memory to preserve multi-systemhomeostasis. <i>J. Clin Invest.</i> Advance online publication doi: 10.1172/JCI121361
ChIP-Seq	BRD4, H3K27Ac	JCI Insight	2018	Anisley Valenciaga et al. Transcriptional targeting of oncogene addiction in medullary thyroid cancer. <i>JCI Insight.</i> 3: e122225
RNA-Seq	NA	PAIN	2018	Pradipta Ray et al. 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> 159:1325
ChIP-Seq	ELL2	The Journal of Immunology	2018	Ashley M. Nelson et al. 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> doi: 10.4049/jimmunol.1800557
ATAC-Seq, ChIP-Seq	H3K36me3, H3K27Ac	PLoS One	2018	Ryan K. Shultzaberger et al. Agnostic detection of genomic alterations by holistic DNA structural interrogation. <i>PLoS ONE</i> 13 (11):e02080554
ChIP-Seq	MTA1	Cancer Medicine	2018	Nasir A. Butt et al. Targeting MTA1/HIF-1 α signaling by pterostilbene in combination with histone deacetylase inhibitor attenuates prostate cancer progression. <i>Cancer Med.</i> 6: 2673
ChIP-Seq	CREB, SWc	Scientific Reports	2018	Jason L. Larabee et al. Unique, Intersecting, and Overlapping Roles of C/EBP β and CREB in Cells of the Innate Immune System <i>Sci Rep.</i> 8:16931
ChIP-Seq, Next-Gen Bisulfite-Seq	H3K4me3, RNA Pol II	Agronomy	2018	Martin Schmidt et al. Methylome and epialleles in rice epilines selected for energy use efficiency. <i>Agronomy.</i> 8: 163
ChIP-Seq	BRD4	Atherosclerosis	2018	Dean Gilham et al. Apabetalone down regulates factors and pathways associated with vascular calcification <i>Atherosclerosis</i> 280:75
ChIP-Seq	GATA2	The American Journal of Human Genetics	2018	Katelyn M. Mika et al. An ancientfecundability-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 et al. 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

Technique	Target	Journal	Year	Reference
MeDIP-Seq	5-Methylcytosine	Scientific Reports	2018	Jose P. Silva et al. 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 et al. Phosphorylation state of ZFP24 controls oligodendrocyte differentiation. <i>Cell Reports.</i> 23: 2254
ChIP-Seq, Histone PTM Quantitation	H3K9me3, H3K27me3	Cell Reports	2018	Jessica Camacho et al. 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 et al. 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 et al. Macrophage/microglial Ezh2 facilitates autoimmune inflammation through inhibition of <i>Socs3</i> . <i>J Exp Med.</i> 215: 1365
ChIP-Seq	STAT3, STAT5, ROCK2	Scientific Reports	2018	Wei Chen et al. 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 et al. Mice lacking the transcriptional regulator <i>Bhlhe40</i> 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 et al. 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 et al. 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 et al. 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 et al. 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 et al. 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 et al. TRRAP is a central regulator of human multiciliated cell formation. <i>J Cell Biol.</i> 217: 1941
ChIP-Seq	BRD4	Blood	2018	Rebecca Kohnken et al. Diminished microRNA-29b level is associated with BRD4 mediated activation of oncogenes in cutaneous T-cell lymphoma. <i>Blood.</i> 131: 771

Technique	Target	Journal	Year	Reference
RNA-Seq	NA	Cell Communication and Signaling	2018	Keri Callegari et al. 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 et al. MTA drives malignant progression and bone metastasis in prostate cancer. <i>Mol. Oncol.</i> 12:1596
ChIP-Seq	FOXG1	Molecular Neurobiology	2018	Stefan Weise et al. FOXG1 Regulates PRKAR2B Transcriptionally and Posttranscriptionally via miR200 in the Adult Hippocampus. <i>Mol. Neurobio.</i> S12035-018-1444.
ChIP-Seq	RNA Pol II, H2A.Z, CDK9, ANP32e	Biochimica et Biophysica Acta	2018	Hyewon Shin et al. Transcriptional regulation mediated by H2A.Z via ANP32e-dependent inhibition of protein phosphatase 2A. <i>Biochim Biophys Acta.</i> 1861: 481
ChIP-Seq	H3K4me3, H3K-9me3, H3K27me3	Proceedings of the Royal Society B	2018	Theresa K. Kelly et al. 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 et al. 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
ChIP-Seq	H3K27Me3	G3 - Genes Genomics Genetics	2018	James Ferguson et al. 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 et al. Origin and differentiation of human memory CD8 T cells after vaccination. <i>Nature.</i> 552: 362
ChIP-Seq	10 different Histone mods, CTCF, RNA Pol II, BRD4	Nature	2017	Elizabeth Stewart et al. Orthotopic patient-derived xenografts of paediatric solid tumours. <i>Nature.</i> 549: 96
ChIP-Seq	BACH2	Nature Communications	2017	Nicolas Hipp et al. 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, PPARα	Cell Reports	2017	Alison Iroz et al. A specific ChREBP and PPARα 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 et al. 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 et al. 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 et al. Defining baseline epigenetic landscapes in the rat liver. <i>Epigenomics.</i> 9: 1503
ChIP-Seq	H3K36me3	The Journal of Clinical Investigation	2017	Huairui Yuan et al. 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 et al. Breast cancer suppression by progesterone receptors is mediated by their modulation of estrogen receptors and RNA polymerase III. <i>Cancer Res.</i> 77: 4934

Technique	Target	Journal	Year	Reference
ChIP-Seq	EZH2, H3K27me3, H3K4me3	Molecular Cancer Therapy	2017	Dorothy Brach et al. 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 et al. 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 et al. 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 et al. 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 et al. Role of ERα 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 et al. Repression of stress-induced LINE-1 expression protects cancer cell subpopulations from lethal drug exposure. <i>Cancer Cell.</i> 32: 1
ChIP-Seq	BRD4, H3K9me3, RNA Pol II	Neuron	2017	Issam Aldiri et al. The dynamic epigenetic landscape of the retina during development, reprogramming and tumorigenesis. <i>Neuron.</i> 94: 550
ChIP-Seq	LMX1b	Development	2017	Endika Haro et al. Lmx1b-targeted cis-regulatory modules involved in limb dorsalization. <i>Dev.</i> 444: 2009
ChIP-Seq	EZH2	Proceedings of the National Academy of Sciences	2017	Yongfeng Liu et al. Epithelial EZH2 serves as an epigenetic determinant in experimental colitis by inhibiting TNFα-mediated inflammation and apoptosis. <i>Proc Natl Acad Sci.</i> 114: E3796
ChIP-Seq	Estrogen Receptor	Endocrinology	2017	Laurel A. Coons et al. DNA sequence constraints define functionally active steroid nuclear receptor binding sites in chromatin. <i>Endocrinology.</i> 158: 3212
ChIP-Seq	ETV5	Proceedings of the National Academy of Sciences	2017	Zhen Zhang et al. 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 et al. 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 data analysis	H3K27me3	Scientific Reports	2017	Brid O'Leary et al. 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 et al. 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 et al. 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 et al. TRIM28 controls a gene regulatory network based on endogenous retroviruses in human neural progenitor cells. <i>Cell Reports</i> 18: 1

Technique	Target	Journal	Year	Reference
ChIP-Seq	CBX3/HP1	Scientific Reports	2017	Michael Sun et al. Cbx3/HP1 γ deficiency confers enhanced tumor-killing capacity on CD8+ T cells. <i>Sci Rep.</i> 7: 42888
ChIP-Seq	H3K27Ac	Scientific Reports	2017	Yao Shen et al. 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 et al. 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 et al. Neuronal activity, TGF β -Signaling and unpredictable chronic stress modulate transcription of Gadd45 family members and DNA methylation in the hippocampus. <i>Cereb Cortex.</i> 27: 4166
MeDIP-Seq	NA	Genes, Brain and Behavior	2017	M. Fonteneau et al. Inhibition of DNA methyltransferases regulates cocaine self-administration by rats: a genome-wide DNA methylation study. <i>Genes Brain Behav.</i> 16:313
ChIP-Seq	SRC-2	PLoS Genetics	2017	Shruthy Suresh et al. SRC-2-mediated coactivation of anti-tumorigenic target genes suppresses MYC-induced liver cancer. <i>PLoS Genet.</i> 13: e1006650
ChIP-Seq	Progesterone Receptor	Biology of Reproduction	2017	Margeaux Wetendorf et al. Decreased epithelial progesterone receptor A at the window of receptivity is required for preparation of the endometrium for embryo attachment. <i>Biol Reprod.</i> 96: 313.
ChIP-Seq	SLY1	Cell Death and Differentiation	2017	Charlotte Moretti et al. SLY regulates genes involved in chromatin remodeling and interacts with TBL1XR1 during sperm differentiation. <i>Cell Death Differ.</i> 24: 1029
ChIP-Seq	G34R	Acta Neuropathologica Communications	2017	Farhana Haque et al. 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 et al. 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 et al. 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 et al. 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 et al. Pan-cancer analysis of somatic copy-number alterations implicates IRS4 and IGF2 in enhancer hijacking. <i>Nat Genet.</i> 49: 65
ChIP-Seq	OTX2, MITF, BRD4, H3K27Ac	Cancer Cell	2016	Pascal D. Johann et al. 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 et al. 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 et al. Non-Cell-Autonomous Regulation of Prostate Epithelial Homeostasis by <i>Androgen Receptor</i> . <i>Mol Cell.</i> 63: 976
ChIP-Seq, Spike-in	BRD4	The Journal of Clinical Investigation	2016	Mark L. McClelland et al. CCAT1 is an enhancer-templated RNA that predicts BET sensitivity in colorectal cancer. <i>J Clin Invest.</i> 126: 639

Technique	Target	Journal	Year	Reference
hMeDIP-Seq	5-Hydroxymethyl-cytosine	Nature Communications	2016	Carolina M. Greco et al. 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 et al. 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 et al. 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 et al. 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 et al. Mouse Dux is myotoxic and shares partial functional homology with its human paralog DUX4. <i>Hum Mol Genet.</i> 25: 4577
ChIP-Seq	H3K27me3	Molecular and Cellular Biology	2016	Kyung Hyun Yoo et al. Histone demethylase KDM6A controls the mammary luminal lineage through enzyme-independent mechanisms. <i>Mol Cell Biol.</i> 36: 2108
ChIP-Seq	H3K27Ac	Nature Genetics	2016	Pratiti Bandopadhyay et al. 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 et al. 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 et al. 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, H3K-27me3, HLX, LHX2, LMX1A	Nature	2016	Charles Y. Lin et al. 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 et al. 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 et al. 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 et al. Loss of Tet1 associated 5-hydroxymethylcytosine is concomitant with aberrant promoter hypermethylation in liver cancer. <i>Cancer Res.</i> 76: 3097
ChIP-Seq, ChIP-qPCR	H3K36me3	Oncogene	2016	Thai H. Ho et al. 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 et al. 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 et al. Antimicrobial peptide LL-37 participates in the transcriptional regulation of melanoma cells. <i>J Cancer.</i> 26: 2341

Technique	Target	Journal	Year	Reference
ChIP-Seq, ChIP-qPCR	p53	Molecular Oncology	2016	Cheryl Chan et al. 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 et al. The regulatory repertoire of PLZF and SALL4 in undifferentiated spermatogonia. <i>Development.</i> 143:1893
ChIP-Seq	H3K27Ac, RNA Pol II, JDP2	Genome Research	2016	Petros Kolovos et al. Binding of nuclear factor κB to noncanonical consensus sites reveals its multimodal role during the early inflammatory response. <i>Genome Res.</i> 26:1478
ChIP-Seq	PLZF	PLoS Genetics	2016	Ramakrishna Kommagani et al. The Promyelocytic leukemia zinc finger transcription factor Is critical for human endometrial stromal cell decidualization. <i>PLoS Genet.</i> 12: e1005937
ChIP libraries, sequencing, analysis	H3K27me3	Developmental Dynamics	2016	Oyvind Dahle et al. 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	Androgen Receptor, Estrogen Receptor	Molecular Cancer Research	2016	Nicholas D'Amato et al. Cooperative Dynamics of AR and ER Activity in Breast Cancer. <i>Mol Cancer Res.</i> 14:1054
ChIP-Seq	NR4A1	PLoS One	2016	Ryan P. Duren et al. 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
ChIP-Seq	MTA1	Oncotarget	2016	Swati Dhar et al. Dietary pterostilbene is a novel MTA1-targeted chemopreventive and therapeutic agent in prostate cancer. <i>Oncotarget.</i> 7: 18469
ChIP-Seq	Myc-tagged PET-1	The Journal of Neuroscience	2016	Steven C. Wyler et al. 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 et al. TGF-β3-expressing CD4 ⁺ CD25 ⁺ LAG3 ⁺ regulatory T cells control humoral immune responses. <i>Nat Commun.</i> 6: 6329
ChIP-Seq	HA-tagged SHOX2	Development	2015	Wenduo Ye et al. 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	BAF180/PBRM1	Molecular Cell	2015	Bokai Zhu et al. 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 et al. 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 et al. 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 et al. 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 Genet.</i> 11: e1005693
ChIP-Seq	Progesterone Receptor	Cell Reports	2015	Vincent J. Lynch et al. Ancient transposable elements transformed the uterine regulatory landscape and transcriptome during the evolution of mammalian pregnancy. <i>Cell Rep.</i> 10: 551

Technique	Target	Journal	Year	Reference
ChIP-Seq	EGR1	PLoS One	2015	Anthony A. Portale et al. 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 et al. 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 et al. 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	NA	The Journal of Clinical Investigation	2015	Coralie Hoareau-Aveilla et al. Reversal of microRNA-150 silencing disadvantages crizotinib-resistant NPM-ALK(+) cell growth. <i>J Clin Invest.</i> 125: 3505
ChIP-Seq, ChIP-qPCR	β -Arrestin-1	Cancer Research	2015	Smitha Pillai et al. β -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, TFIIB	Circulation: Heart Failure	2015	Danish Sayed et al. Acute targeting of general transcription factor IIB restricts cardiac hypertrophy via selective inhibition of gene transcription. <i>Circ Heart Fail.</i> 8: 138
ChIP-Seq	H3K4me3	Journal of the National Cancer Institute	2015	Cristian Taccioli et al. 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 et al. 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 et al. Lhx6 and Lhx8 promote palate development through negative regulation of a cell cycle inhibitor gene, p57Kip2. <i>Hum Mol Genet.</i> 24: 5024
ChIP-Seq	SOX9	Nucleic Acids Research	2015	Zhongcheng Shi et al. Context-specific role of SOX9 in NF- κ B mediated gene regulation in colorectal cancer cells. <i>Nucleic Acids Res.</i> 43: 6257
ChIP-Seq	H3K27Me3	Nucleic Acids Research	2015	Kyung Hyun Yoo et al. 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 et al. 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 et al. 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 et al. 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 et al. Progesterone receptor transcriptome and cistrome in decidualized human endometrial stromal cells. <i>Endocrinology.</i> 156: 2239
ChIP-Seq	H3K27me3, H3K4me1, H3K4me3, H3K27Ac, p300, RAR α	Immunity	2015	Chrysothemis C. Brown et al. Retinoic acid is essential for Th1 cell lineage stability and prevents transition to a Th17 cell program. <i>Immunity.</i> 42: 1

Technique	Target	Journal	Year	Reference
ChIP-Seq	SIX1	Cancer Cell	2015	Jenny Wegert et al. 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 et al. 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 et al. Specific phosphorylation of histone demethylase KDM3A determines target gene expression in response to heat shock. <i>PLOS Biol</i> . 12: e1002026
ChIP-Seq	SRC-1	Molecular Endocrinology	2014	Mounia Tannour-Louet et al. 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 et al. 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> . 15: 761
ChIP-Seq	GATA2, FOXA1, p300, CBP, SRC-1, SRC-2, SRC-3, Androgen Receptor	Proceedings of the National Academy of Sciences	2014	Bin He et al. GATA2 facilitates steroid receptor coactivator recruitment to the androgen receptor complex. <i>Proc Natl Acad Sci</i> . 111: 18261
Next-Gen Bisulfite-Seq	NA	Cell Reports	2014	Deepak Kumar et al. Fibroblast growth factor maintains chondrogenic potential of limb bud mesenchymal cells by modulating DNMT3A recruitment. <i>Cell Rep</i> . 8: 1419
ChIP-Seq	PPAR γ	Biology of Reproduction	2014	Kelsey E. Brooks et al. Peroxisome proliferator activator receptor gamma (PPAR γ) regulates conceptus elongation in sheep. <i>Biol Reprod</i> . 92: 42.
ChIP-Seq, LightSwitch	PXR, p300, H3K4me1, H3K27Ac	PLoS Genetics	2014	Robin P. Smith et al. 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 et al. 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	SRC-3	The Journal of Clinical Investigation	2014	Jun Qin et al. 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 et al. Identification of a face enhancer reveals direct regulation of LIM homeobox 8 (Lhx8) by wingless-int (WNT)/ β -catenin Signaling. <i>J Biol Chem</i> . 289: 30289
ChIP-Seq, ChIP-qPCR	ZBTB20	Journal of Medical Genetics	2014	Malene B. Rasmussen et al. 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 et al. Enhancer hijacking activates GF11 family oncogenes in medulloblastoma. <i>Nature</i> . 511: 428
ChIP-Seq	H3K4me3, H3K27me3	Nature	2014	Volker Hovestadt et al. Decoding the regulatory landscape of medulloblastoma using DNA methylation sequencing. <i>Nature</i> . 510: 537

Technique	Target	Journal	Year	Reference
ChIP-qPCR	EZH2, SUZ12, H3K27me3	Molecular Cancer Therapy	2014	Sarah K. Knutson et al. 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 α , RAR β , RXR α	American Journal of Physiology - Gastrointestinal and Liver Physiology	2014	Yuqi He et al. 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 et al. 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 et al. Novel DNA motif binding activity observed In vivo with an estrogen receptor a mutant mouse. <i>Mol Endocrinol.</i> 26: 899
ChIP-Seq	H3K27me3	Bioinformatics	2014	Yanxiao Zhang et al. 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 et al. 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 et al. SRC-2 Is an essential coactivator for orchestrating metabolism and circadian rhythm. <i>Cell Rep.</i> 6: 633
Next-Gen Bisulfite-Seq	NA	Nature Neuroscience	2014	Brian G. Dias et al. 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 et al. 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 et al. 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 et al. PARP-14 binds specific DNA sequences to promote Th2 cell gene expression. <i>PLoS One.</i> 8: e83127
Next-Gen Bisulfite-Seq	NA	Clinical Cancer Research	2013	David S. Shames et al. 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	CCND1 (Cyclin D1)	Cancer Research	2013	Xiaoming Ju et al. 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 et al. 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 et al. 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 et al. COUP-TFII regulates human endometrial stromal genes involved in inflammation. <i>Mol Endocrinol.</i> 27: 2041

Technique	Target	Journal	Year	Reference
ChIP-Seq	GFI1	Nature Immunology	2013	Chauncey J. Spooner et al. 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 et al. PARP-14 and its enzyme activity regulates Th2 differentiation and allergic airway disease. <i>J Allergy Clin Immunol.</i> 131: 521
ChIP-Seq	Progesterone Receptor	Molecular Endocrinology	2013	Ashlee R. Lain et al. Research Resource: Progesterone receptor targetome underlying mammary gland branching morphogenesis. <i>Mol Endocrine.</i> 27: 1743
ChIP-Seq	PPAR α	Chemico-Biological Interactions	2013	Patrick D. McMullen et al. A map of the PPAR α transcription regulatory network for primary human hepatocytes. <i>Chem Biol Interact.</i> 209: 14
ChIP-Seq	H3K36me3	Cancer Discovery	2013	Lynn Bjerke et al. 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 et al. 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 et al. 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 et al. 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 et al. 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 α , RNA Pol II	PLoS One	2013	Astrid Kosters et al. Sexually dimorphic genome-wide binding of retinoid X receptor alpha (RXR α) determines male-female differences in the expression of hepatic lipid processing genes in mice. <i>PLoS One.</i> 8: e71538
ChIP-qPCR	NF κ B (p52 and p65)	Journal of Biological Chemistry	2013	Sarah L. Doyle et al. Nuclear factor κ B2 p52 protein has a role in antiviral immunity through I κ B kinase epsilon-dependent induction of Sp1 protein and interleukin 15. <i>J Biol Chem.</i> 288: 25066
ChIP-chip, ChIP-qPCR	C/EBP β , RNA Pol II	Journal of Biological Chemistry	2013	Hana Vakili et al. CCAAT-enhancer-binding protein β (C/EBP β) 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 et al. 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 et al. Vitamin D receptor ChIP-seq in primary CD4+ 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 et al. Atrial identity is determined by a COUP-TFII regulatory network. <i>Dev Cell.</i> 25: 417
ChIP-Seq	H3K4me3	PLoS One	2013	Alex Gutteridge et al. 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 et al. 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

Technique	Target	Journal	Year	Reference
ChIP-Seq	LSD1, H3K4me2	Molecular and Cellular Biology	2012	Venugopalan D. Nair et al. 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 et al. EZH2 inhibition as a therapeutic strategy for lymphoma with EZH2-activating mutations. <i>Nature.</i> 492: 108
ChIP-Seq data analysis	pTyr37 H2B	Nature Structural and Molecular Biology	2012	Kiran Mahajan et al. H2B Tyr37 phosphorylation suppresses expression of replication-dependent core histone genes. <i>Nat Struct Mol Biol.</i> 19: 930
ChIP-chip	B-MYB	PLoS One	2012	Ming Zhan et al. 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-tagged BAP1, OGT, HCF1	Science	2012	Anwasha Dey et al. 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 et al. 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 et al. 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 et al. 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 et al. Whole-Genome Estrogen Receptor Binding in Mouse Uterine Tissue Revealed by ChIP-Seq. <i>Mol Endocrinol.</i> 26: 887
ChIP-qPCR	PDX1, TCF3	Chemistry & Biology	2012	Alice Kiselyuk et al. 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 et al. 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 et al. 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 et al. GATA4 expression is primarily regulated via a miR-26b-dependent post-transcriptional mechanism during cardiac hypertrophy. <i>Cardiovasc Res.</i> 93: 645
ChIP-Seq	Farnesoid X Receptor	American Journal of Physiology - Gastrointestinal and Liver Physiology	2012	Julia Yue Cui et al. 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 et al. 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-tagged CCND1 (Cyclin D1)	The Journal of Clinical Investigation	2012	Mathew C. Casimiro et al. ChIP sequencing of cyclin D1 reveals a transcriptional role in chromosomal instability in mice. <i>J Clin Invest.</i> 122: 833

Technique	Target	Journal	Year	Reference
ChIP-Seq	RBPJ	Stem Cells	2012	Yaochen Li et al. 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
ChIP-chip	Progesterone Receptor	The FASEB Journal	2011	Heather L. Franco et al. Epithelial progesterone receptor exhibits pleiotropic roles in uterine development and function. <i>FASEB J</i> . 26: 1218
ChIP-Seq	MSGN1	Nature Communications	2011	Ravindra B Chalamalasetty et al. The Wnt3a/ β -catenin target gene Mesogenin1 controls the segmentation clock by activating a Notch signalling program. <i>Nat Commun</i> . 2: 390
ChIP-qPCR	IRF4, cMyc	British Journal of Haematology	2011	Antonia Lopez-Girona et al. (2011) 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 et al. Increased expression of IRF4 and ETS1 in CD4+ cells from patients with intermittent allergic rhinitis. <i>Allergy</i> . 67: 33
ChIP-Seq	NELF-b (COBRA1)	Journal of Biological Chemistry	2011	Jianlong Sun et al. 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 et al. 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 et al. Targeting MYC dependence in cancer by inhibiting BET bromodomains. <i>Proc Natl Acad Sci</i> . 108:16669
ChIP-chip	C/EBP β	Journal of Molecular Endocrinology	2011	Aristides Lytras et al. 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 Yudt et al. 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 et al. The Transcription Factor 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 et al. 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 et al. 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
Next-Gen Bisulfite-Seq	NA	PLoS One	2011	Michela Deleidi et al. 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	BMC Biology	2011	Genevieve Lavoie et al. PKC isoforms interact with and phosphorylate DNMT1. <i>BMC Biol</i> . 9: 31
ChIP-qPCR	Aryl Hydrocarbon Receptor	Toxicology and Applied Pharmacology	2011	K Nadira De Abrew et al. 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
ChIP-qPCR	PPAR γ	Investigative Ophthalmology and Visual Science	2011	Gerard A Rodrigues et al. Differential Effects of PPAR γ Ligands on Oxidative Stress-Induced Death of Retinal Pigmented Epithelial Cells. <i>Invest Ophthalmol Vis Sci</i> . 52: 890

Technique	Target	Journal	Year	Reference
ChIP-chip	p73	Proceedings of the National Academy of Sciences	2011	Jennifer M Rosenbluth et al. 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 et al. 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	Aryl Hydrocarbon Receptor, cMaf	Nature Immunology	2010	Lionel Apetoh et al. 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 et al. 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 et al. 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 et al. 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	Pregnane X Receptor	Toxicological Sciences	2010	Julia Yue Cui et al. Genetic and Epigenetic Regulation and Expression Signatures of Glutathione S-Transferases in Developing Mouse Liver. <i>Toxicol Sci.</i> 116: 32
ChIP-Seq	SRC-3, RNA Pol II	Molecular Endocrinology	2010	Rainer B Lanz et al. 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 et al. 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> 286: 15380
ChIP-chip	Aryl Hydrocarbon Receptor	Toxicological Sciences	2010	K Nadira De Abrewet al. 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 et al. 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 et al. The Androgen Receptor Modulates Expression of Genes with Critical Roles in Muscle Development and Function. <i>Mol Endocrinol.</i> 24: 1665
ChIP-qPCR	NFkB (p50 and p65)	Cardiovascular Research	2010	Alina G Sofronescu et al. FGF-16 is a target for adrenergic stimulation through NF-kB activation in postnatal cardiac cells and adult mouse heart. <i>Cardiovasc Res.</i> 87: 102
ChIP-Seq	Pregnane X Receptor	Nucleic Acids Research	2010	Julia Yue Cui et al. 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 et al. Temporal ChIP-on-Chip of RNA-Polymerase-II to detect novel gene activation events during photoreceptor maturation. <i>Mol Vis.</i> 16: 252
ChIP-chip	STAT4	The Journal of Immunology	2009	Seth R Good et al. 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 et al. A Gene Signature-Based Approach Identifies mTOR as a Regulator of p73. <i>Mol Cell Biol.</i> 28: 5951
ChIP-qPCR	Aryl Hydrocarbon Receptor	Nature	2008	Francisco J Quintana et al. Control of Treg and TH17 cell differentiation by the aryl hydrocarbon receptor. <i>Nature.</i> 453:65