

**Chromatin Modifiers**  
**DNA Methylation**  
**Histones & Oncohistones**  
**Transcription Regulation**

## Highly Characterized Antibodies for Epigenetics and Gene Regulation

At Active Motif, we are committed to providing the highest-quality antibodies for studying epigenetics in the context of histone and DNA modifications. Our antibodies are manufactured in-house, and undergo rigorous development and validation procedures to ensure their quality and performance. Our extensive, novel portfolio of monoclonal, polyclonal and AbFlex® recombinant antibodies are validated for the applications you need – chromatin immunoprecipitation (ChIP), ChIP-Seq, CUT&Tag, Western blot, and more.

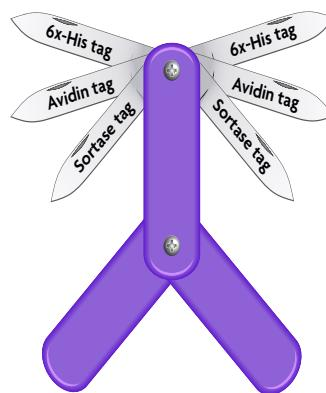
To browse our complete list of antibodies and application data, please visit [activemotif.com/abs](http://activemotif.com/abs)

### AbFlex® Recombinant Antibodies

#### Uniquely designed for multiple labeling methods

AbFlex® antibodies are recombinant antibodies (rAbs) that have been generated using defined DNA sequences to produce highly specific, reproducible antibodies. The unique advantages of the AbFlex® antibody are its flexible labeling and purification options. Each AbFlex® antibody contains a Sortase recognition motif (LPXTG) to covalently add fluorophores, enzymatic substrates (e.g., HRP), peptides, DNA, drugs, or other labels to the antibody in a directed and reproducible manner. Every antibody also contains a 6xHis tag, which can be used with nickel-based purification systems, and an avidin tag sequence for enzymatic biotin conjugation using the biotin ligase, BirA.

**AbFlex is available  
for a range of  
Histone Markers,  
DNA Methylation and  
Transcription Factors**

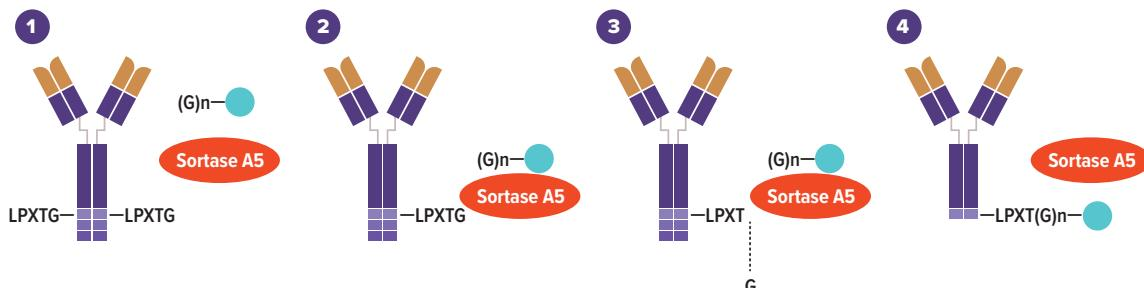


**AbFlex™**

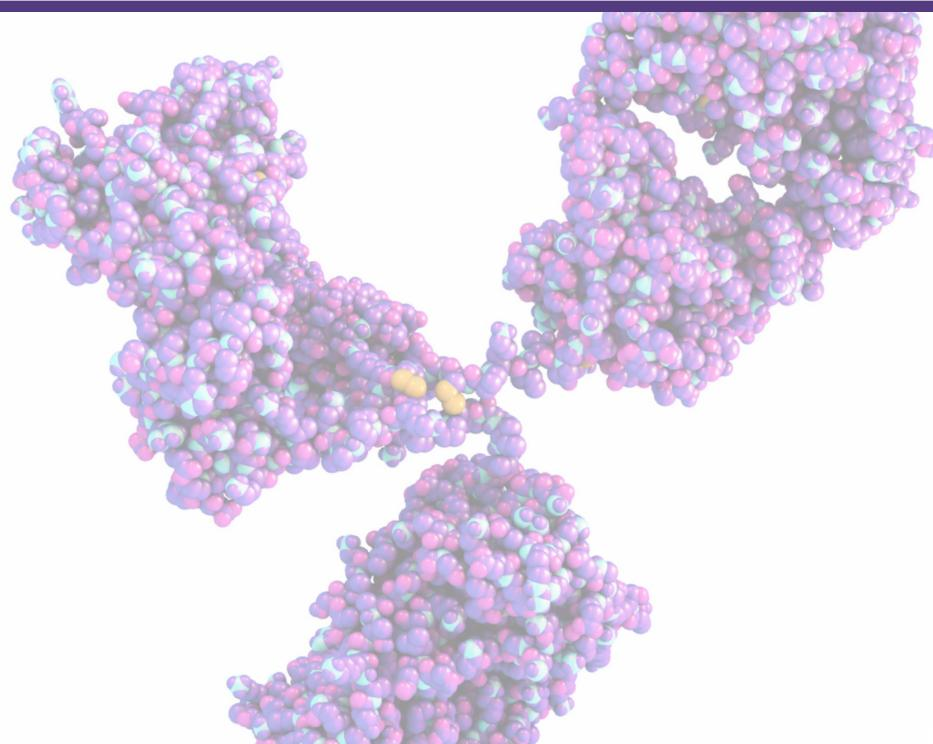
## Sortag-IT™ Labeling Kits

### For site specific conjugation to AbFlex® rAbs

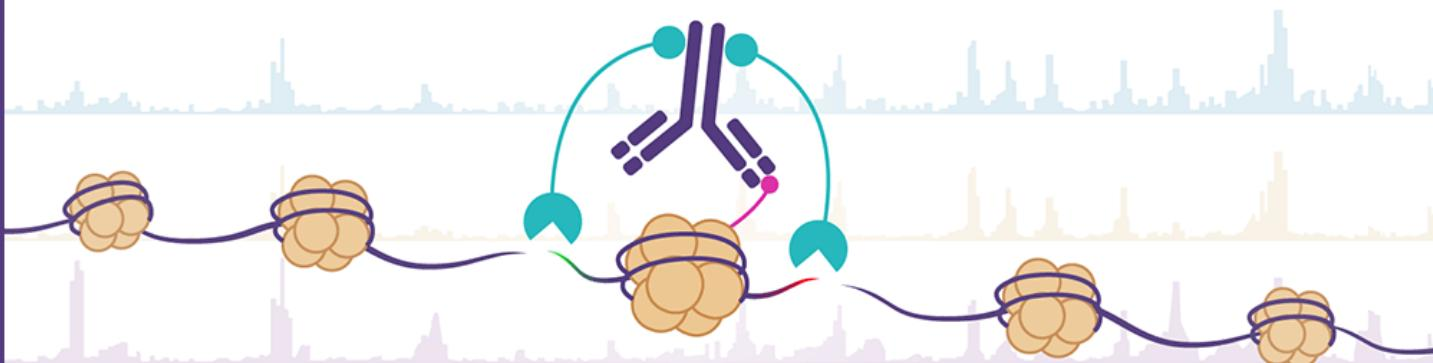
Sortag-IT™ Labeling Kits are designed for site-specific labeling of Active Motif's highly specific AbFlex® recombinant antibodies (rAb) via the Sortase tag recognition sequence (LPXTG) that is incorporated into the heavy chains of each AbFlex antibody. Directly conjugate fluorophores, enzymes (HRP, AP), biotin, peptides, DNA, carbohydrates or other labels to the terminal end of the heavy chain to ensure the antigen binding site remains available for target recognition.

**Figure:**

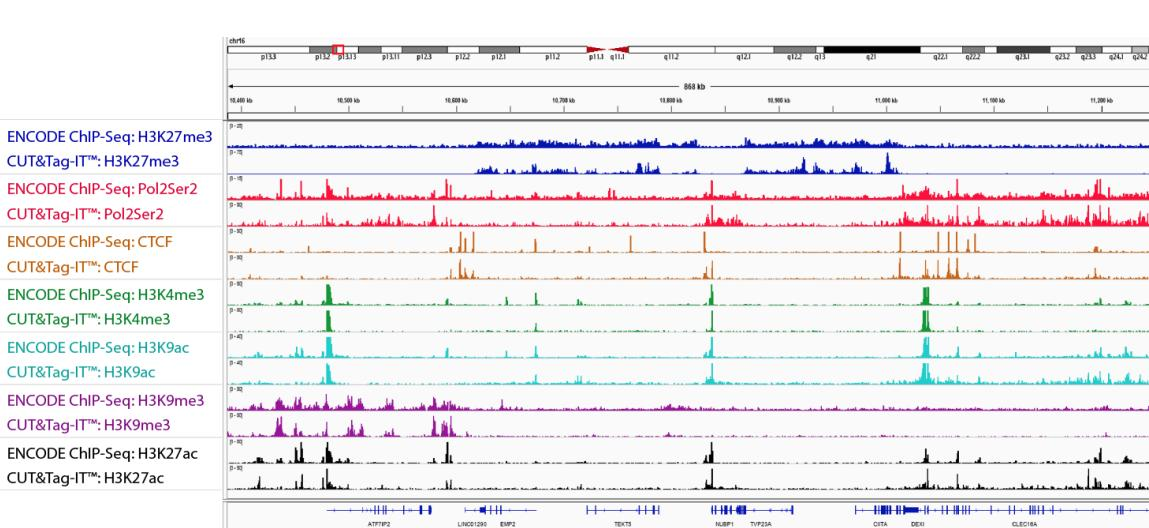
Sortase A belongs to a family of transpeptidases found in Gram-positive bacteria that catalyze the attachment of poly-Glycine (G)n labels to the sortase recognition sequence. The Sortag-IT Labeling Kits use Active Motif's Sortase A5 pentamutant from *Staphylococcus aureus* that is significantly more active than wild-type Sortase to provide a faster, more efficient labeling reaction. Simply add your AbFlex antibody with the poly-Glycine label, add Sortase A5 and incubate for approximately 1 hour. Purification columns are included to remove excess label and Stop Solution is provided to inactivate the Sortase A5 enzyme. Labeled antibodies are ready for downstream analysis.



## CUT&Tag-Validated Antibodies



**Cleavage Under Target and Tagmentation (CUT&Tag)** is an alternative method to traditional ChIP assays that can reveal interactions between histone modifications and other proteins and the DNA they interact with or to identify DNA binding sites for proteins of interest. The CUT&Tag assay requires a highly specific primary antibody to the protein target of interest that must specifically detect the protein in its native, unfixed form. Active Motif specializes in manufacturing high-quality antibodies to histones, histone modifications, chromatin proteins, and other factors, including a growing list of antibodies that we have experimentally validated in-house to work well in CUT&Tag assays.



**Figure: CUT&Tag-IT Assay Kit data correlate well with ENCODE ChIP-Seq data.**

Chromatin landscapes from human K562 cells are shown across a 868 kb segment of the human genome region (Chr16: 10,400-11,250) for H3K27me3 (dark blue), phospho-Pol2Ser2 (red), CTCF (orange), H3K4me3 (green), H3K9ac (light blue) H3K9me3 (purple), and H3K27ac (black) showing high concordance of CUT&Tag-IT Assay Kit signals from 100,000 cells compared with ENCODE ChIP-Seq data from at least one million cells. Antibodies used in the CUT&Tag assays for the following targets were: H3K27me3: Cat. No 39155, CTCF: Cat. No. 61311, H3K4me3: Cat. No. 39159, H3K9ac: Cat. No. 39917, H3K9me3: Cat. No. 39161, H3K27ac: Cat. No. 39133.

Description	Applications	Cat. No.	Description	Applications	Cat. No.
3-Methylcytosine (3-mC) antibody (pAb)	DB, ICC, IF	61111	beta Tubulin antibody (mAb)	IHC, WB	61157
3-Methylcytosine (3-mC) antibody (pAb)	DB, ICC, IF	61179	beta-Catenin antibody (pAb)	IP, WB	61487
5-Carboxylcytosine (5-caC) antibody (pAb)	DB, ICC, IF, IHC, WB	61225	BLIMP1 / PRDM1 antibody (pAb)	IF, WB	61053
5-Carboxylcytosine (5-caC) antibody (pAb)	DB, ICC, IF	61229	BMI-1 antibody (mAb)	ChIP, ChIP-Seq, IF, IHC, WB	39993
5-Formylcytosine (5-fC) antibody (pAb)	DB, ICC, IF, IHC, WB	61223	BMI-1 antibody (mAb)	ICC, IF, WB	61089
5-Formylcytosine (5-fC) antibody (pAb)	DB, ICC, IF, WB	61227	AbFlex® BMI-1 antibody (rAb)	ChIP, ChIP-Seq, WB	91195
5-Hydroxymethylcytosine (5-hmC) antibody (mAb)	DB, ICC, IF, IHC, MeDIP	39999	Boris / CTCFL antibody (pAb)	ChIP, ChIP-Seq, WB	39851
5-Hydroxymethylcytosine (5-hmC) antibody (pAb)	DB, FC, ICC, IF, IHC, MeDIP	39769	BRD2 antibody (pAb)	ChIP, ChIP-Seq	61797
5-Hydroxymethylcytosine (5-hmC) antibody (pAb)	DB, IF, IHC, MeDIP	39791	BRD3 antibody (pAb)	ChIP, ChIP-Seq, WB	61489
5-mCpA antibody (mAb)	DB	61783	AbFlex® BRD4 antibody (rAb)	ChIP, ChIP-Seq	91301
5-Methylcytosine (5-mC) antibody (mAb)	DB, IF, MeDIP	61479	BRD8 / SMAP2 antibody (pAb)	ChIP, ChIP-Seq, WB	61007
AbFlex® 5-methylcytosine antibody (rAb)	DB	91187	BRD9 antibody (pAb)	IP, WB	65705
AATF / Che-1 antibody (mAb)	IF, WB	61789	BrdU antibody (mAb)	IHC	61273
Acetyl-Lysine antibody (rAb)	ICC, WB	91315	AbFlex® BrdU antibody (rAb)	IF	91269
ACTL6B antibody (pAb)	WB	61691	Bridging Antibody for Mouse IgG	ChIP, IP	53017
AbFlex® Ago 1/2/3 antibody (rAb)	IP	91217	BRPF1 antibody (pAb)	WB	61541
Ago1 antibody (mAb)	IP, WB	61071	Bub1 antibody (mAb)	ICC, IF	39735
Ago1/2/3 antibody (mAb)	ChIP, ICC, IF, IHC, IP, WB	39937	BubR1 antibody (mAb)	ICC, IF	39747
Ago2 antibody (mAb)	ICC, IF, IP, WB	39853	c-Fos antibody (pAb)	WB	61421
Ago3 antibody (mAb)	ChIP, ICC, IF, IP	39787	c-Jun antibody (pAb)	WB	65707
Ago4 antibody (mAb)	IP, WB	39855	c-Myc antibody (mAb)	EMSA, IP, WB	61075
Ago4 antibody (pAb)	WB	39813	AbFlex® c-Myc antibody (rAb)	WB	91213
AIB1 / SRC-3 antibody (pAb)	ChIP, WB	39797	C17orf96 antibody (pAb)	ChIP-Seq, WB	61753
AIB3 antibody (pAb)	IHC, WB	39835	C17orf96 phospho Ser15 antibody (pAb)	DB	61745
AID antibody (mAb)	ChIP, WB	39885	CABIN1 antibody (pAb)	IP, WB	61597
Aiolos antibody (mAb)	ChIP, WB	39657	CAF-1 p150 antibody (mAb)	ICC, IF, IP, WB	39997
alpha Tubulin antibody (mAb)	ICC, IF, WB	39527	CAF-1 p60 antibody (mAb)	ICC, IF, IP, WB	39995
AbFlex® AM-Tag antibody (rAb)	ChIP, ELISA, WB	91111	CARM1 antibody (pAb)	ChIP, IHC, IP, WB	39251
Androgen Receptor antibody (pAb)	ChIP, ChIP-Seq, WB	39781	Cas9 antibody (mAb)	ChIP, ICC, IF, IP, WB	61577
ARID3A antibody (pAb)	WB	61615	Cas9 antibody (mAb)	ChIP, IF, IP, WB	61757
ASH2L antibody (mAb)	ICC, IF, IHC, IP, WB	61359	AbFlex® Cas9 antibody (rAb)	ChIP, ChIP-Seq, ELISA, WB	91123
ASH2L antibody (pAb)	ChIP, ChIP-Seq, WB	39099	Caspase-3 antibody (mAb)	WB	40924
ATAD2 antibody (pAb)	IP, WB	61369	CBX8 antibody (pAb)	WB	61237
ATF1 antibody (pAb)	WB	61563	CDK8 antibody (pAb)	ChIP, ChIP-Seq, WB	61481
ATF4 antibody (pAb)	WB	61655	CDK9 antibody (pAb)	WB	61559
ATM phospho Ser1981 antibody (mAb)	ICC, IF, WB	39529	CDKN2A antibody (pAb)	IP, WB	61619
AbFlex® ATM phospho Ser1981 antibody (rAb)	DB, WB	91207	CDKN2A p14ARF antibody (pAb)	WB	61689
ATRX antibody (pAb)	WB	61049	CENP-A phospho Ser18 antibody (pAb)	DB, ICC, IF, WB	61483
Aurora A antibody (mAb)	FC, ICC, IF, IP, WB	61709	CENP-B antibody (pAb)	ICC, IF	61183
Aurora B antibody (pAb)	ICC, IF, WB	39261	CENP-B antibody (pAb)	ICC, IF, IP, WB	61287
BAP1 antibody (pAb)	WB	61501	CENP-E antibody (mAb)	ICC, IF	39619
BAZ2A antibody (pAb)	IF, WB	61707	CHD1 antibody (mAb)	ICC, IF, WB	39989
BCL11A antibody (mAb)	ICC, IF, IHC, WB	61231	CHD1 antibody (pAb)	WB	39729
Bcl6 antibody (mAb)	IHC, WB	61193	CHD2 antibody (pAb)	WB	39363
AbFlex® Bcl6 antibody (rAb)	WB	91287	CHD3 antibody (mAb)	ICC, IF, IP, WB	61469
Bcl7A antibody (pAb)	IP, WB	65695	AbFlex® CHD3 antibody (rAb)	IP, WB	91273

The lists above are a small subset of our antibodies, we also offer each in a **convenient sample size**.

For an up-to-date list of all available antibodies, please visit [activemotif.com/abs](http://activemotif.com/abs).

Description	Applications	Cat. No.	Description	Applications	Cat. No.
Chd5 antibody (mAb)	ICC, IF, WB	61683	FEN1 antibody (pAb)	WB	65709
CID antibody (mAb)	ICC, IF	61735	FosB antibody (mAb)	WB	40960
CoREST antibody (pAb)	WB	39955	FOXG1 antibody (pAb)	ChIP, ChIP-Seq, WB	61211
CTBP antibody (pAb)	WB	39847	FOXL2 antibody (pAb)	WB	61283
CTBP2 antibody (pAb)	ChIP, ChIP-Seq, WB	61261	FOXM1 antibody (pAb)	WB	61693
CTCF antibody (pAb)	ChIP, ChIP-Seq, CUT&Tag, EMSA, ICC, IF, IHC, WB	61311	FOXO1/FKHR antibody (pAb)	IP, WB	39629
AbFlex® CTCF antibody (rAb)	ChIP-Seq, CUT&Tag, WB	91285	FOXP1 antibody (mAb)	ChIP, EMSA, WB	61309
CtlP antibody (mAb)	ChIP, ICC, IF, IP, WB	61141	FTO antibody (pAb)	WB	61431
CUL4B antibody (pAb)	WB	61445	GATA-4 antibody (pAb)	WB	39893
Cxxc4 antibody (pAb)	WB	61679	GATA-6 antibody (pAb)	ICC, IF, WB	61063
DAP-3 antibody (mAb)	WB	40963	Geminin antibody (pAb)	IF, IP, WB	61135
DAX-1 / NR0B1 antibody (mAb)	ICC, IF, IHC, WB	39983	GFI1 antibody (pAb)	WB	61553
DDB1 antibody (pAb)	WB	61127	GLI1 antibody (pAb)	WB	61215
Ddit3 antibody (mAb)	IP, WB	61349	GLIS1 antibody (mAb)	IF, WB	61801
DHX9 antibody (mAb)	ICC, IF, WB	61705	GLYR1 antibody (pAb)	WB	65693
DMAP1 antibody (pAb)	IP, WB	61675	Goosecoid antibody (pAb)	WB	61121
DNA-RNA Hybrid antibody (mAb)	ChIP, ChIP-Seq, DB, EMSA, ICC, IF, IP	65683	HA-Tag antibody (mAb)	IP	39627
DNMT1 antibody (mAb)	ChIP, IHC, IP, WB	39204	AbFlex® HA-Tag antibody (rAb)	IP	91271
DNMT1 antibody (mAb)	ChIP, IP, WB	61467	HDAC1 antibody (mAb)	WB	39881
DNMT1 antibody (pAb)	IF, IP, WB	39905	HDAC1 antibody (mAb)	ChIP, ICC, IF, IHC, IP, WB	39531
AbFlex® DNMT1 antibody (rAb)	WB	91175	HDAC1 antibody (pAb)	ChIP, ChIP-Seq, WB	40967
DNMT1 monomethyl Lys142 antibody (pAb)	WB	61497	AbFlex® HDAC1 antibody (rAb)	ChIP-Seq, WB	91215
DNMT3A antibody (mAb)	ChIP, IHC, WB	39206	HDAC11 antibody (pAb)	WB	39208
DNMT3B antibody (mAb)	WB	61503	HDAC2 antibody (mAb)	ChIP, ICC, IF, WB	39533
DNMT3B antibody (pAb)	WB	39899	AbFlex® HDAC2 antibody (rAb)	ChIP-Seq, WB	91197
DOT1L antibody (pAb)	WB	39953	HDAC3 antibody (mAb)	IP, WB	39717
Drosha antibody (pAb)	IP, WB	39783	HDAC3 antibody (pAb)	ChIP, WB	40968
EED antibody (mAb)	ChIP, ChIP-Seq, IHC, WB	61203	AbFlex® HDAC3 antibody (rAb)	WB	91199
AbFlex® EED antibody (rAb)	ChIP, ChIP-Seq, ELISA	91135	HDAC4 antibody (pAb)	ChIP, ChIP-Seq, WB	40969
AbFlex® EHMT2 (G9a) antibody (rAb)	WB	91229	HDAC5 antibody (pAb)	ChIP, IP, WB	40970
EKLF antibody (mAb)	ChIP, IP, WB	61233	HDAC6 antibody (pAb)	ChIP, WB	40971
Elk-1 antibody (pAb)	IP, WB	61301	HDAC9 antibody (pAb)	WB	39793
ELP3 antibody (pAb)	WB	39949	HIF-1 alpha antibody (mAb)	IHC, WB	61275
ERR-α antibody (pAb)	WB	61047	HIF-1 alpha antibody (pAb)	ChIP, ChIP-Seq, WB	39665
Estrogen Receptor α antibody (mAb)	ChIP, IP, WB	61035	AbFlex® HIF-1 alpha antibody (rAb)	ELISA	91139
Estrogen Receptor α antibody (pAb)	ChIP, WB	65715	HIPK2 antibody (mAb)	ICC, IF, IHC, IP, WB	39677
AbFlex® Estrogen Receptor α antibody (rAb)	WB	91411	HIRA antibody (mAb)	ICC, IF, WB	39557
Estrogen Receptor β antibody (mAb)	IHC, IP, WB	39767	HIRA antibody (mAb)	IP, WB	39555
ETO / RUNX1T1 antibody (pAb)	IP, WB	61109	AbFlex® His-Tag antibody (rAb)	WB	91183
Ets-1 antibody (pAb)	ChIP-Seq, WB	39580	Histone H1 antibody (mAb)	ChIP, ChIP-Seq, IF, IP, WB	61785
EuHMT1 antibody (pAb)	WB	39777	Histone H1 antibody (pAb)	ICC, IF, WB	39707
EZH1 antibody (pAb)	WB	61583	Histone H1 antibody (pAb)	WB	61201
EZH2 antibody (mAb)	ChIP	39875	Histone H1.0 antibody (pAb)	WB	61417
EZH2 antibody (pAb)	ChIP, ChIP-Seq, IP	39901	Histone H1.5S17ph antibody (pAb)	DB, WB	61107
EZH2 antibody (pAb)	ChIP	61641	Histone H2A antibody (pAb)	WB	39209
AbFlex® EZH2 antibody (rAb)	IP	91159	Histone H2A antibody (pAb)	ChIP, IP, WB	39235

The lists above are a small subset of our antibodies, we also offer each in a **convenient sample size**.

For an up-to-date list of all available antibodies, please visit [activemotif.com/abs](http://activemotif.com/abs).

Description	Applications	Cat. No.	Description	Applications	Cat. No.
Histone H2A antibody (pAb)	ChIP, WB	39945	Histone H3.3S31ph antibody (pAb)	DB, IF, WB	39637
AbFlex® Histone H2A antibody (rAb)	WB	91325	Histone H3.cs1 antibody (pAb)	WB	39573
Histone H2A, acidic patch antibody (pAb)	ChIP, WB	39111	Histone H3.X/Y antibody (mAb)	ChIP, ICC, IF, WB	61161
Histone H2A, C-terminal antibody (pAb)	IF, WB	39591	Histone H3ac (pan-acetyl) antibody (pAb)	ChIP, ChIP-Seq, DB, WB	39139
Histone H2A.J antibody (pAb)	DB, WB	61793	Histone H3ac (pan-acetyl) antibody (pAb)	ChIP, DB, IF, WB	61637
Histone H2A.X antibody (pAb)	ChIP, ICC, IF, WB	39689	Histone H3K122me1 antibody (pAb)	DB, WB	39367
Histone H2A.XS139ph antibody (pAb)	DB, ICC, IF, WB	39117	Histone H3K14ac antibody (mAb)	ChIP, ChIP-Seq, WB	61433
AbFlex® Histone H2A.XS139ph antibody (rAb)	WB	91395	Histone H3K14ac antibody (pAb)	ChIP, DB, WB	39697
Histone H2A.XS139ph antibody (rAb)	ICC, WB	91313	Histone H3K14ac antibody (pAb)	ChIP, DB, WB	65701
Histone H2A.XY142ph antibody (pAb)	DB, WB	61617	AbFlex® Histone H3K14ac antibody (rAb)	ELISA, WB	91397
Histone H2A.Z antibody (pAb)	ChIP, ChIP-Seq, DB, IF, WB	39113	Histone H3K14bz antibody (pAb)	DB, WB	39303
Histone H2A.Z antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39943	Histone H3K14me1 antibody (pAb)	DB, IF, WB	61509
Histone H2A.Z.2.2 antibody (mAb)	WB	61737	Histone H3K14me2 antibody (pAb)	WB	39349
Histone H2A/H4S1ph antibody (pAb)	DB, WB	39115	Histone H3K14me3 antibody (pAb)	DB, WB	61549
Histone H2AK119ub antibody (pAb)	DB, WB	65697	Histone H3K18ac antibody (pAb)	ChIP, DB, ICC, IF, WB	39129
Histone H2AK5ac antibody (pAb)	ChIP, DB, IF, WB	39107	Histone H3K18ac antibody (pAb)	ChIP, DB, ICC, IF, IHC, WB	39587
Histone H2AK9ac antibody (pAb)	DB, IF, WB	39109	Histone H3K18ac antibody (pAb)	ChIP, DB, ICC, IF, WB	39693
Histone H2AQ104me1 antibody (pAb)	DB, WB	61605	Histone H3K18ac antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39755
Histone H2AS129ph antibody (pAb)	ChIP, DB, WB	39271	Histone H3K18me1 antibody (pAb)	DB, WB	39667
Histone H2AT120ph antibody (pAb)	DB, WB	39391	Histone H3K23ac antibody (pAb)	ChIP, DB, ICC, IF, IP, WB	39131
Histone H2AT120ph antibody (pAb)	DB, IF, WB	61195	Histone H3K23me1 antibody (pAb)	DB, ICC, IF, WB	39387
Histone H2Av antibody (mAb)	ChIP, WB	61751	Histone H3K23me2 antibody (pAb)	ChIP, DB, ICC, IF, WB	39653
Histone H2B antibody (pAb)	WB	39210	Histone H3K23me3 antibody (pAb)	DB, ICC, IF	61499
Histone H2B antibody (pAb)	WB	39125	Histone H3K23pr antibody (pAb)	DB, WB	61397
Histone H2B antibody (pAb)	ChIP, WB	39237	Histone H3K27ac antibody (mAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39685
Histone H2B antibody (pAb)	WB	39947	Histone H3K27ac antibody (pAb)	ChIP, ChIP-Seq, CUT&Tag, DB, ICC, IF, IHC, WB	39133
Histone H2BK120ac antibody (pAb)	ChIP, ChIP-Seq, DB, IF, WB	39119	AbFlex® Histone H3K27ac antibody (rAb)	ChIP-Seq, CUT&Tag, ELISA, WB	91193
Histone H2BK120ub1 antibody (mAb)	ChIP, WB	39623	Histone H3K27me1 antibody (mAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	61015
Histone H2BK12ac antibody (pAb)	ChIP, DB, IF, WB	39669	Histone H3K27me1 antibody (pAb)	DB, ICC, IF, WB	39377
Histone H2BK15ac antibody (pAb)	WB	61321	Histone H3K27me1 antibody (pAb)	DB, WB	39889
Histone H2BK16ac antibody (pAb)	ChIP, DB, IF, WB	39121	Histone H3K27me2 antibody (mAb)	ChIP, ChIP-Seq, WB	61435
Histone H2BK46ac antibody (pAb)	ChIP, WB	39571	Histone H3K27me2 antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39245
Histone H2BK46me2 antibody (pAb)	DB, ICC, IF, WB	39567	Histone H3K27me2 antibody (pAb)	DB, ICC, IF, WB	39919
Histone H2BK5ac antibody (pAb)	ChIP, ChIP-Seq, DB, IF, WB	39123	Histone H3K27me3 antibody (mAb)	ChIP, ChIP-Seq, IF, WB	39535
Histone H2BK5me1 antibody (pAb)	DB, WB	61439	Histone H3K27me2me3 antibody (mAb)	ChIP, ICC, IF, WB	39536
Histone H2BS14ph antibody (mAb)	ICC, IF, WB	61011	Histone H3K27me3 antibody (mAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	61017
Histone H3 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, WB	39763	Histone H3K27me3 antibody (pAb)	ChIP, ChIP-Seq, CUT&Tag, DB, ICC, IF, IHC, WB	39155
Histone H3 antibody (mAb)	ICC, IF, WB	61647	Histone H3K27me3 antibody (pAb)	ChIP, CUT&Tag, DB, ICC, IF, IP, WB	39156
Histone H3 antibody (pAb)	ChIP-Seq, WB	61799	AbFlex® Histone H3K27me3 antibody (rAb)	DB, WB	91167
Histone H3, C-terminal antibody (pAb)	ELISA, WB	39451	AbFlex® Histone H3K27me3 antibody (rAb)	ChIP, DB, WB	91403
AbFlex® Histone H3, C-terminal antibody (rAb, Mouse)	WB	91297	Histone H3K36ac antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39379
AbFlex® Histone H3, C-terminal antibody (rAb, Rabbit)	ChIP, WB	91299	Histone H3K36me0 antibody (pAb)	DB, WB	65111
Histone H3.1 / 3.2 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, WB	61629	Histone H3K36me1 antibody (pAb)	DB, WB	61351
Histone H3.1S28ph antibody (mAb)	ICC, IF, WB	61697	Histone H3K36me2 antibody (mAb)	ChIP, DB, IF, WB	61019
AbFlex® Histone H3.3 antibody (rAb)	ChIP-Seq, WB	91191	Histone H3K36me2 antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39255
Histone H3.3K27M antibody (pAb)	WB	61803			

The lists above are a small subset of our antibodies, we also offer each in a **convenient sample size**.

For an up-to-date list of all available antibodies, please visit [activemotif.com/abs](http://activemotif.com/abs).

Description	Applications	Cat. No.
Histone H3K36me2 antibody (pAb)	DB, WB	39891
Histone H3K36me3 antibody (mAb)	ChIP, ChIP-Seq, DB, IF, WB	61021
Histone H3K36me3 antibody (pAb)	ChIP, ChIP-Seq, CUT&Tag, DB, WB	61101
AbFlex® Histone H3K36me3 antibody (rAb)	ChIP, ChIP-Seq, WB	91265
Histone H3K37ac antibody (pAb)	DB, IF, WB	61587
Histone H3K4ac antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39381
Histone H3K4ac antibody (pAb)	ChIP, ChIP-Seq, DB, WB	61809
Histone H3K4K9K18 (pan-biotinylated) antibody (pAb)	DB, ICC, IF, WB	61391
AbFlex® Histone H3K4me0 antibody (rAb, mouse)	DB, WB	91267
AbFlex® Histone H3K4me0 antibody (rAb, rabbit)	DB, WB	91317
Histone H3K4me1 antibody (mAb)	ChIP, ChIP-Seq, DB, WB	39635
Histone H3K4me1 antibody (pAb)	ChIP, DB, WB	61633
Histone H3K4me1 antibody (pAb)	ChIP, ChIP-Seq, WB	39299
Histone H3K4me1 antibody (pAb)	ChIP, ChIP-Seq, WB	61781
AbFlex® Histone H3K4me1 antibody (rAb)	WB	91289
Histone H3K4me2 antibody (mAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39679
Histone H3K4me2 antibody (pAb)	ChIP, ChIP-Seq, CUT&Tag, DB, WB	39141
Histone H3K4me2 antibody (pAb)	ChIP, DB, ICC, IF, WB	39913
AbFlex® Histone H3K4me2 antibody (rAb)	ChIP, ChIP-Seq, WB	91321
Histone H3K4me3 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, WB	61379
Histone H3K4me3 antibody (pAb)	ChIP, ChIP-Seq, CUT&Tag, DB, ICC, IF, IHC, WB	39159
Histone H3K4me3 antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39915
AbFlex® Histone H3K4me3 antibody (rAb)	ChIP, ChIP-Seq, DB, WB	91263
Histone H3K56ac antibody (mAb)	ChIP, DB	61061
Histone H3K56ac antibody (pAb)	ChIP, ChIP-Seq, DB, WB	39281
AbFlex® Histone H3K56ac antibody (rAb)	ChIP-Seq, ELISA	91127
Histone H3K56me1 antibody (pAb)	DB, WB	39273
Histone H3K56me2 antibody (pAb)	ChIP, DB, WB	39277
Histone H3K79ac antibody (pAb)	ChIP, DB, WB	39565
Histone H3K79me1 antibody (pAb)	ChIP, DB, WB	39145
Histone H3K79me1 antibody (pAb)	ChIP, DB, WB	39921
AbFlex® Histone H3K79me2 antibody (rAb)	ChIP, DB, WB	91399
Histone H3K79me2 antibody (pAb)	ChIP, ChIP-Seq, DB, WB	39143
Histone H3K79me2 antibody (pAb)	ChIP, DB, WB	39923
Histone H3K9ac antibody (mAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	61251
Histone H3K9ac antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, WB	61663
Histone H3K9ac antibody (pAb)	ChIP, ChIP-Seq, DB, IF, WB	39137
Histone H3K9ac antibody (pAb)	DB, ICC, IF, WB	39585
Histone H3K9ac antibody (pAb)	ChIP, ChIP-Seq, CUT&Tag, DB, ICC, IF, WB	39917
AbFlex® Histone H3K9ac antibody (rAb)	ChIP, ChIP-Seq, ELISA, ICC, IF, WB	91103
Histone H3K9me0 antibody (mAb)	WB	61399
AbFlex® Histone H3K9me0 antibody (rAb)	ChIP, ChIP-Seq, ELISA, ICC, IF, WB	91155
Histone H3K9me1 antibody (mAb)	ChIP, DB, WB	39681
Histone H3K9me1 antibody (pAb)	ChIP, ChIP-Seq, DB, WB	39887
Histone H3K9me1 antibody (pAb)	ChIP-Seq, DB, ICC, IF, WB	39249
Histone H3K9me1me2me3 antibody (pAb)	DB, ICC, IF, WB	39241

Description	Applications	Cat. No.
Histone H3K9me2 antibody (mAb)	ChIP, DB, ICC, IF, WB	39683
Histone H3K9me2 antibody (pAb)	ChIP, DB, ICC, IF, WB	39239
Histone H3K9me2 antibody (pAb)	ChIP, DB, ICC, IF, WB	39375
Histone H3K9me2 antibody (pAb)	ChIP, DB, ICC, IF, WB	39753
AbFlex® Histone H3K9me2 antibody (rAb)	WB	91335
Histone H3K9me3 antibody (mAb)	ChIP, DB, ICC, IF, WB	61013
Histone H3K9me3 antibody (mAb)	DB, ICC, IF, WB	39285
Histone H3K9me3 antibody (pAb)	ChIP, ChIP-Seq, CUT&Tag, DB, ICC, IF, IHC, WB	39161
Histone H3K9me3 antibody (pAb)	ChIP, CUT&Tag, DB, ICC, IF, WB	39765
Histone H3R17me2a (asymmetric) antibody (pAb)	ChIP, DB, IF, IHC, WB	39709
Histone H3R17me2a18ac antibody (pAb)	ChIP, DB, WB	61613
Histone H3R8me1 antibody (pAb)	DB, WB	39673
Histone H3R8me2a (asymmetric) antibody (pAb)	ChIP, ChIP-Seq, DB, WB	39651
Histone H3S10ph antibody (mAb)	ELISA, ICC, IF, WB	61623
Histone H3S10ph antibody (mAb)	DB, ICC, IF, WB	39636
Histone H3S10ph antibody (pAb)	ChIP, DB, ICC, IF, WB	39253
AbFlex® Histone H3S10ph antibody (rAb)	DB, WB	91131
Histone H3S10phS28ph antibody (pAb)	ChIP, ICC, IF, WB	39147
Histone H3S28ph antibody (mAb)	ICC, IF, WB	61695
Histone H3S28ph antibody (mAb)	IF, IHC	39098
Histone H3S28ph antibody (mAb)	WB	61325
Histone H3T11ph antibody (pAb)	ChIP, ChIP-Seq, DB, WB	39151
Histone H3T11ph antibody (pAb)	DB, WB	61197
Histone H3T32ph antibody (mAb)	ICC, IF, WB	61777
Histone H3T3ph antibody (pAb)	DB, ICC, IF, WB	39153
Histone H3T45ph antibody (pAb)	DB, WB	39737
Histone H4 antibody (mAb)	ChIP, WB	61521
Histone H4 antibody (pAb)	ChIP, WB	39269
Histone H4 antibody (pAb)	WB	61199
Histone H4 antibody (pAb)	ChIP, ChIP-Seq, WB	61299
AbFlex® Histone H4 antibody (rAb)	WB	91295
Histone H4ac (pan-acetyl) antibody (mAb)	ICC, IF, IHC, WB	39967
Histone H4ac (pan-acetyl) antibody (pAb)	ChIP, DB, ICC, IF, WB	39243
Histone H4ac (pan-acetyl) antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, IHC, WB	39925
Histone H4H18ph antibody (pAb)	DB, WB	61411
Histone H4H18ph antibody (pAb)	DB, WB	61413
Histone H4K12ac antibody (mAb)	ChIP, DB, WB	61527
Histone H4K12ac antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39165
Histone H4K12ac antibody (pAb)	ChIP, DB, ICC, IF, WB	39927
Histone H4K16ac antibody (mAb)	ChIP, DB, IF, WB	61529
Histone H4K16ac antibody (pAb)	ChIP, ChIP-Seq, DB, IF, WB	39167
Histone H4K16ac antibody (pAb)	ChIP, DB, IHC, WB	39929
Histone H4K20ac antibody (mAb)	ChIP, DB	61531
Histone H4K20me1 antibody (mAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	39727
AbFlex® Histone H4K20me1 antibody (rAb)	WB	91291
Histone H4K20me2 antibody (mAb)	WB	39539

The lists above are a small subset of our antibodies, we also offer each in a **convenient sample size**.

For an up-to-date list of all available antibodies, please visit [activemotif.com/abs](http://activemotif.com/abs).

Description	Applications	Cat. No.
Histone H4K20me2 antibody (mAb)	DB, WB	61533
Histone H4K20me3 antibody (mAb)	ChIP, DB, ICC, WB	39671
AbFlex® Histone H4K20me3 antibody (rAb)	ChIP, ChIP-Seq, ELISA, ICC, IF, WB	91107
Histone H4K31me1 antibody (pAb)	DB, IF, WB	39385
Histone H4K5ac antibody (mAb)	ChIP, DB, ICC, IF, WB	61523
Histone H4K5ac antibody (pAb)	ChIP, DB, ICC, IF, WB	39169
Histone H4K5ac antibody (pAb)	ChIP, DB, ICC, IF, WB	39583
Histone H4K5ac antibody (pAb)	ChIP, ChIP-Seq, DB, IF, WB	39699
AbFlex® Histone H4K5ac/K8ac antibody (rAb)	ChIP, ChIP-Seq, DB, ELISA, WB	91319
Histone H4K8ac antibody (mAb)	ChIP, DB, WB	61525
Histone H4K8ac antibody (pAb)	WB	39171
Histone H4K8ac antibody (pAb)	ChIP, ChIP-Seq, DB, ICC, IF, WB	61103
Histone H4K8K12K16 (pan-biotinylated) antibody (pAb)	DB	61393
Histone H4R3me2a (asymmetric) antibody (pAb)	ChIP, DB, IF, WB	39705
Histone H4R3me2s (symmetric) antibody (pAb)	ChIP, ChIP-Seq, DB, WB	61187
Histone macroH2A1 antibody (pAb)	ICC, IF, WB	39593
Histone macroH2A1.1 antibody (pAb)	ICC, IF, IHC, WB	39871
Histone macroH2A1.2 antibody (mAb)	ICC, IF, IHC, WB	61427
Histone macroH2A2 antibody (pAb)	ICC, IF, IHC, WB	39873
HMGA1 antibody (pAb)	ICC, IF, WB	39615
HMGA2 / HMGI-C antibody (pAb)	IP, WB	61041
HMGN1 antibody (pAb)	IP, WB	39833
Hmgn2 antibody (pAb)	WB	61353
HNF-3α / FOXA1 antibody (mAb)	ChIP, ChIP-Seq, IHC, WB	39837
HNF-3β / FOXA2 antibody (pAb)	IHC, WB	39827
HNF4A antibody (pAb)	ICC, IF, WB	61189
HOXA9 antibody (pAb)	WB	39825
HP1 gamma phospho Ser93 antibody (pAb)	WB	39655
HP1α antibody (pAb)	WB	39295
HP1β antibody (mAb)	ICC, IF, IP, WB	39979
HRP Goat anti-Mouse IgG	ELISA, IHC, WB	15014
HRP Goat anti-Rabbit IgG	ELISA, IHC, WB	15015
HSC70 / HSP7C antibody (mAb)	ICC, IF, IP, WB	61207
Htz1 / Histone H2A.Z antibody (pAb)	ChIP, ChIP-Seq, WB	39647
IFNA1 antibody (mAb)	ELISA	61603
Ikaros antibody (mAb)	ChIP, ChIP-Seq, WB	39355
Ikaros antibody (pAb)	ChIP, EMSA, WB	39291
IKK $\alpha$ /IKK $\beta$ antibody (pAb)	WB	40957
IKK $\gamma$ antibody (mAb)	WB	40908
INCENP antibody (mAb)	ICC, IF	39259
ING5 antibody (mAb)	IP, WB	65685
AbFlex® ING5 antibody (rAb)	WB	91329
IRF-2 antibody (pAb)	WB	61281
IRF-3 antibody (pAb)	ChIP, WB	39371
IRF-4 antibody (pAb)	WB	61511
IRF-5 antibody (pAb)	ChIP, ChIP-Seq, WB	61269

Description	Applications	Cat. No.
IκB $\alpha$ antibody (mAb)	WB	40903
IκB $\alpha$ phospho Ser32,36 antibody (mAb)	WB	40904
JMJD2A antibody (mAb)	WB	39815
AbFlex® JMJD2A antibody (rAb)	ELISA, WB	91143
JMJD2B / KDM4B antibody (pAb)	ICC, IF, IP, WB	61221
AbFlex® JMJD2B / KDM4B antibody (rAb)	IP, WB	91305
AbFlex® JMJD2C / KDM4C antibody (rAb)	WB	91307
AbFlex® JMJD2D / KDM4D antibody (rAb)	DB, IP, WB	91369
JMJD2D antibody (pAb)	WB	39247
JMJD3 / KDM6B antibody (mAb)	WB	61387
JMJD6 antibody (pAb)	WB	61493
JunB antibody (pAb)	WB	39549
JunD antibody (pAb)	ChIP-Seq, IP, WB	61403
KAT2A (GCN5) antibody (pAb)	IP, WB	39831
KAT5 antibody (pAb)	WB	61545
KDM1B / LSD2 antibody (pAb)	WB	61457
AbFlex® KDM1B / LSD2 antibody (rAb)	WB	91327
KDM2B antibody (mAb)	WB	65699
KDM4C antibody (pAb)	WB	61687
AbFlex® KDM5A antibody (rAb)	ChIP-Seq, WB	91211
Ki-67 antibody (pAb)	WB	39799
KLF4 antibody (pAb)	WB	39745
KLF5 antibody (pAb)	ChIP, ChIP-Seq, ICC, IF, WB	61099
KLF6 antibody (mAb)	IHC, WB	61297
KLF6 antibody (pAb)	WB	61243
KMT2E (MLL5) antibody (pAb)	WB	61447
L3MBTL1 antibody (pAb)	ChIP, WB	39182
L3MBTL2 antibody (pAb)	IP, WB	39569
Lamin A, mature antibody (mAb)	ICC, IF, IP, WB	39961
Lamin A/C antibody (mAb)	ChIP, ICC, IF, IP, WB	39287
AbFlex® Lamin A/C antibody (rAb)	IF, WB	91189
LAP2 alpha antibody (mAb)	WB	39267
LexA DNA-binding Domain antibody (pAb)	ChIP, WB	39184
LIN28A antibody (pAb)	ICC, IF, WB	61191
LSD1 / KDM1A antibody (pAb)	ChIP, IP, WB	39186
LXR- $\alpha$ antibody (pAb)	ChIP, ChIP-Seq, WB	61175
LXR- $\beta$ antibody (pAb)	ChIP, ChIP-Seq, WB	61177
MAD1 antibody (mAb)	ICC, IF	39659
MAD2L1 antibody (mAb)	WB	39399
MAF antibody (pAb)	WB	61635
MALT1 antibody (mAb)	WB	39395
MASH1 antibody (mAb)	IHC, WB	61271
MAZ antibody (mAb)	ChIP, EMSA, WB	39935
AbFlex® MAZ antibody (rAb)	WB	91205
MBD1 antibody (mAb)	WB	39215
MBD1 antibody (pAb)	WB	39857

The lists above are a small subset of our antibodies, we also offer each in a **convenient sample size**.  
 For an up-to-date list of all available antibodies, please visit [activemotif.com/abs](http://activemotif.com/abs).

Description	Applications	Cat. No.	Description	Applications	Cat. No.
MBD2 antibody (pAb)	WB	39547	NFATC2 antibody (pAb)	IP, WB	61621
MBD3 antibody (mAb)	WB	39216	NFkB p100 antibody (pAb)	WB	39687
MCM2 antibody (pAb)	WB	61717	NFkB p65 antibody (mAb)	ChIP, WB	40916
MeCP2 antibody (mAb)	ChIP, ICC, IF, IHC, IP, WB	61291	NFkB p65 antibody (pAb)	WB	39283
MeCP2 phospho Ser80 antibody (pAb)	WB	39733	NFkB p65 antibody (pAb)	ChIP, ChIP-Seq, EMSA, WB	39369
MED1 antibody (pAb)	IP, WB	61065	NFkB p65 phospho Ser529 antibody (pAb)	WB	39691
MED15 antibody (pAb)	IP, WB	61589	NFkB p65 phospho Ser536 antibody (pAb)	WB	39675
MEIS 1/2/3 antibody (mAb)	ICC, IF	39795	NKK2.5 antibody (pAb)	WB	61267
AbFlex® MEIS 1/2/3 antibody (rAb)	WB	91259	NonO antibody (pAb)	IP, WB	61247
Menin antibody (pAb)	ChIP, ChIP-Seq, WB	61005	Notch1 antibody (mAb)	WB	61147
MEP50 antibody (pAb)	WB	39190	AbFlex® Notch1 antibody (rAb)	WB	91243
METTL1 / WDR4 antibody (pAb)	WB	65113	NR3C1 antibody (pAb)	WB	61519
METTL3 antibody (mAb)	IP, WB	65687	NRF1 antibody (pAb)	WB	61673
AbFlex® METTL3 antibody (rAb)		91333	NRF2 antibody (pAb)	ChIP, ChIP-Seq, WB	61599
MGMT antibody (pAb)	IP, WB	39951	Nucleolin antibody (mAb)	ICC, IF, WB	39541
Mi-2 antibody (mAb)	IF, IP, WB	61463	AbFlex® Nucleolin antibody (rAb)	WB	91231
AbFlex® Mi-2 beta (CHD4) antibody (rAb)	WB	91275	Nucleophosmin antibody (pAb)	WB	39849
Mi-2 beta antibody (mAb)	WB	39695	O-GlcNAc antibody (mAb)	IP, WB	61453
Mili / PiwiL2 antibody (mAb)	ICC, IF, IP, WB	61143	Oct-4 antibody (pAb)	ChIP, ChIP-Seq, ICC, IF, WB	39811
AbFlex® Mili / PiwiL2 antibody (rAb)	WB	91253	OGA/O-GlcNAcase antibody (pAb)	IP, WB	61425
MINA antibody (pAb)	WB	61713	OGT/O-GlcNAc transferase antibody (pAb)	ChIP, ChIP-Seq, WB	61355
MITF antibody (mAb)	ChIP, ChIP-Seq, EMSA, ICC, IF, IP, WB	39789	Osteoprotegerin (OPG) antibody (mAb)	WB	40938
AbFlex® MITF antibody (rAb)	ChIP-Seq	91201	p300 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, IP, WB	61401
MLL / HRX antibody (mAb)	ChIP, IP, WB	39829	p53 antibody (mAb)	ChIP, IP, WB	39553
MLL / HRX antibody (pAb)	ChIP, ChIP-Seq, ICC, IF, IP, WB	61295	p53 antibody (pAb)	WB	61657
MMSET / WHSC1 antibody (mAb)	ICC, IF, WB	39879	AbFlex® p53 antibody (rAb)	WB	91247
MOF / MYST1 antibody (pAb)	ChIP, ICC, IF, WB	61245	PAF1 antibody (pAb)	WB	61567
MRG15 antibody (pAb)	WB	39361	PARP-1 C-terminal antibody (pAb)	ChIP, ChIP-Seq, WB	39561
MTA2 antibody (pAb)	WB	39359	PARP-1 N-terminal antibody (pAb)	ChIP, WB	39559
Myc-Tag antibody (mAb)	WB	39279	PARP-1 N-terminal antibody (pAb)	ChIP-Seq, CUT&Tag, IP, WB	61639
AbFlex® Myc-Tag antibody (rAb)	WB	91203	PARP-2 antibody (pAb)	ChIP, IF, IP, WB	39743
Myf-5 antibody (mAb)	ICC, IF, WB	39801	PAX5 antibody (pAb)	WB	61565
MyoD antibody (mAb)	ICC, IF, WB	39991	PAX6 antibody (pAb)	WB	61611
N-Myc antibody (pAb)	ChIP-Seq, WB	61185	PAX7 antibody (mAb)	ICC, IF, WB	39803
N6-Methyladenosine (m6A) antibody (mAb)	DB, IP	61755	PAX9 antibody (mAb)	IHC, WB	61077
N6-Methyladenosine (m6A) antibody (pAb)	DB	61495	PBRM1 antibody (pAb)	ChIP-Seq, WB	61381
AbFlex® N6-Methyladenosine (m6A) antibody (rAb)	DB, IP	91261	PBX1b antibody (mAb)	ChIP, IHC, WB	61165
Nanog antibody (mAb)	ICC, IF, WB	61627	PCIF1/CAPAM antibody (pAb)	IP, WB	65711
Nanog antibody (pAb)	ChIP, ChIP-Seq, ICC, IF, WB	61419	PCL2 antibody (mAb)	WB	61153
AbFlex® Nanog antibody (rAb)	WB	91251	PDX1 antibody (pAb)	WB	61289
Nap1 antibody (pAb)	IF, WB	39577	PGC1-β antibody (pAb)	WB	39859
NAT10 antibody (pAb)	WB	65703	Phc1 antibody (mAb)	ICC, IF, IP, WB	39723
NCAPH2 antibody (mAb)	ICC, IF, WB	61681	PHF20 antibody (pAb)	WB	61123
NCOA1 antibody (pAb)	WB	61653	PHF6 antibody (pAb)	WB	61807
NCoA4 / ELE1 antibody (pAb)	WB	39895	PHF8 antibody (pAb)	WB	39711
NF-1A antibody (pAb)	ICC, IF, IHC, WB	39397	Pht1 / Histone H2A.Z antibody (pAb)	WB	39640

The lists above are a small subset of our antibodies, we also offer each in a **convenient sample size**.

For an up-to-date list of all available antibodies, please visit [activemotif.com/abs](http://activemotif.com/abs).

Description	Applications	Cat. No.	Description	Applications	Cat. No.
Ph1 / Histone H2A.Zac (pan-acetyl) antibody (pAb)	DB, WB	39642	RNA pol II CTD phospho Ser7 antibody (mAb)	ICC, IF, WB	61703
PIM2 antibody (mAb)	IHC, WB	61159	RNA pol II CTD phospho Thr4 antibody (mAb)	ChIP, DB, ELISA, ICC, IF, IP, WB	61361
AbFlex® Pim2 antibody (rAb)	WB	91239	RNA pol II CTD phospho Thr4 antibody (pAb)	WB	61307
PITX2 antibody (pAb)	WB	61303	RNA pol II CTD phospho Tyr1 antibody (mAb)	ChIP, ICC, IF, IP, WB	61383
PKC-B antibody (pAb)	WB	61733	AbFlex® RNA Pol II CTD phospho Tyr1 antibody (rAb)	ChIP-Seq, ELISA, IF, WB	91219
PLZF antibody (mAb)	ChIP, ICC, IF, IP, WB	39987	RNA pol II CTD Ser2ph / Ser5ph antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, WB	61665
POLR2B antibody (pAb)	IP, WB	61557	RXR-γ antibody (mAb)	EMSA, IHC, WB	61033
PP2A antibody (pAb)	ChIP, WB	39192	RYBP antibody (pAb)	WB	61461
PPM1D antibody (pAb)	WB	61591	SaCas9 antibody (mAb)	WB	61787
PRDM1 / BLIMP1 antibody (mAb)	IHC, IP, WB	61167	AbFlex® SaCas9 antibody (rAb)	WB	91279
PRMT2 antibody (pAb)	ICC, IF, WB	61491	SALL4 antibody (pAb)	IP, WB	39957
PRMT5 antibody (pAb)	ChIP-Seq, WB	61001	SAP30 antibody (pAb)	ChIP, ChIP-Seq, WB	39731
PRMT6 antibody (pAb)	IHC, WB	61003	SARS-CoV-2 Spike Antibody (AM001414)	ELISA, NEU	91361
Progerin antibody (mAb)	ICC, IF, WB	39965	SARS-CoV-2 Spike Antibody (AM002413)	ELISA	91379
Prox1 antibody (mAb)	IF, WB	61091	SARS-CoV-2 Spike Antibody (AM002414)	ELISA, NEU	91349
Prox1 antibody (mAb)	ICC, IF, IHC	61093	SARS-CoV-2 Spike Antibody (AM004414)	ELISA, WB	91345
PXR antibody (pAb)	WB	39589	SARS-CoV-2 Spike Antibody (AM005415)	ELISA	91347
Rad21 antibody (mAb)	ICC, IF, WB	39383	SARS-CoV-2 Spike Antibody (AM005505)	ELISA	91383
AbFlex® Rad21 antibody (rAb)	IF, WB	91245	SARS-CoV-2 Spike Antibody (AM006415)	ELISA	91351
Rad51 antibody (pAb)	ICC, IF	39194	SARS-CoV-2 Spike Antibody (AM009105)	ELISA	91337
RAR-α antibody (mAb)	EMSA, IHC, WB	39971	SARS-CoV-2 Spike Antibody (AM015553)	ELISA, NEU	91377
RAR-β antibody (mAb)	EMSA, IHC, WB	61027	SARS-CoV-2 Spike Antibody (AM017553)	ELISA	91387
RAR-γ antibody (mAb)	EMSA, IHC, WB	39973	SARS-CoV-2 Spike Antibody (AM018553)	ELISA	91389
RB1 antibody (pAb)	IP, WB	61585	SARS-CoV-2 Spike Antibody (AM020553)	ELISA	91381
RbAp46/48 antibody (pAb)	ChIP, WB	39198	SARS-CoV-2 Spike Antibody (AM027553)	ELISA	91391
RBBP5 antibody (pAb)	ChIP, ChIP-Seq, WB	61405	SARS-CoV-2 Spike Antibody (AM032414)	ELISA	91359
RBM39 antibody (pAb)	WB	61669	SARS-CoV-2 Spike Antibody (AM038105)	ELISA	91339
RBPJ antibody (mAb)	WB	61073	SARS-CoV-2 Spike Antibody (AM043105)	ELISA	91341
RBPJ antibody (mAb)	ChIP, EMSA, ICC, IF	61505	SARS-CoV-2 Spike Antibody (AM063553)	ELISA	91393
RING1 antibody (pAb)	WB	61517	AbFlex® SARS-CoV-2 Spike Antibody (rAb) (AM001414)	ELISA, NEU	91375
Ring1B antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, WB	39663	AbFlex® SARS-CoV-2 Spike Antibody (rAb) (AM002414)	ELISA, NEU	91363
RNA pol II antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, WB	39097	AbFlex® SARS-CoV-2 Spike Antibody (rAb) (AM004414)	ELISA	91373
RNA pol II antibody (mAb)	ChIP, IF, IP, WB	61081	AbFlex® SARS-CoV-2 Spike Antibody (rAb) (AM009105)	ELISA	91365
RNA pol II antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, WB	61667	AbFlex® SARS-CoV-2 Spike Antibody (rAb) (AM038105)	ELISA	91371
AbFlex® RNA Pol II antibody (rAb)	ChIP, ChIP-Seq, IF, WB	91151	AbFlex® SARS-CoV-2 Spike Antibody (rAb) (AM043105)	ELISA	91367
RNA pol II CTD monomethyl Lys7 antibody (mAb)	ChIP, ChIP-Seq, IP, WB	65691	SARS-CoV-2 Spike IgG/IgM Antibody (AM043105)	ELISA	91385
RNA pol II CTD phospho Ser2 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, IP, WB	61083	SATB1 antibody (pAb)	WB	39839
RNA pol II CTD phospho Ser2 antibody (mAb)	ICC, IF, WB	61699	SATB2 antibody (pAb)	WB	61575
RNA pol II CTD phospho Ser2 antibody (pAb)	IP, WB	39563	SET8 / PR-SET7 antibody (pAb)	WB	61009
AbFlex® RNA pol II CTD phospho Ser2 antibody (rAb)	ChIP, ChIP-Seq, ELISA, ICC, IF, WB	91115	SETD7 / SET7 antibody (pAb)	WB	61451
RNA pol II CTD phospho Ser5 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, IP, WB	61085	SFMBT1 antibody (pAb)	ChIP, ChIP-Seq, IF, IP, WB	61759
RNA pol II CTD phospho Ser5 antibody (mAb)	ICC, IF, WB	61701	SIN3A antibody (pAb)	ChIP, WB	39865
RNA pol II CTD phospho Ser5 antibody (pAb)	ChIP, ICC, IF, WB	39233	SIP1 antibody (mAb)	ICC, IF, IHC, WB	61095
RNA pol II CTD phospho Ser5 antibody (pAb)	ChIP, ChIP-Seq, WB	39749	SIP1 antibody (pAb)	WB	61113
AbFlex® RNA pol II CTD phospho Ser5 antibody (rAb)	ChIP, ChIP-Seq, ELISA, ICC, IF, WB	91119	SIRT1 antibody (mAb)	ICC, IF, IP, WB	39353
RNA pol II CTD phospho Ser7 antibody (mAb)	ChIP, ChIP-Seq, ICC, IF, IP, WB	61087	SIRT2 antibody (pAb)	ICC, IF, WB	61045

The lists above are a small subset of our antibodies, we also offer each in a **convenient sample size**.

For an up-to-date list of all available antibodies, please visit [activemotif.com/abs](http://activemotif.com/abs).

Description	Applications	Cat. No.	Description	Applications	Cat. No.
SIRT2 antibody (pAb)	ICC, IF, WB	61043	Supt5H/SPT5 antibody (mAb)	ChIP-Seq, IP, WB	65689
SIRT2 antibody (pAb)	IP, WB	39903	AbFlex® Supt5H/SPT5 antibody (rAb)	IP, WB	91331
SIRT2 phospho Ser331 antibody (mAb)	WB	61363	SUV39H1 antibody (mAb)	ChIP, ICC, IF, IP, WB	39785
SIRT6 antibody (pAb)	IP, WB	39911	Suz12 antibody (mAb)	ChIP, IHC, IP	39877
SMAD3 antibody (pAb)	WB	61249	Suz12 antibody (pAb)	ChIP, ChIP-Seq, IF, WB	39357
SMAD4 antibody (pAb)	IP, WB	61631	AbFlex® Suz12 antibody (rAb)	IP	91163
SMARCA1 / SNF2L1 antibody (mAb)	IHC, WB	61465	TAZ / WWTR1 antibody (pAb)	WB	61265
AbFlex® SMARCA1 / SNF2L1 antibody (rAb)	WB	91225	TBK1 antibody (mAb)	WB	40909
SMARCA2 / BRM antibody (mAb)	ChIP, ICC, IF, WB	39805	TCF4 antibody (pAb)	WB	61609
SMARCA4 / BRG1 antibody (mAb)	ICC, IF, WB	39807	TCF7 antibody (pAb)	WB	61579
SMARCA4 antibody (mAb)	ICC, IF, WB	61625	TCF7L1 / TCF3 antibody (pAb)	ChIP, ChIP-Seq, WB	61125
SMARCAL1 / HARP antibody (mAb)	IP, WB	61115	TDG / Thymine-DNA glycosylase antibody (pAb)	WB	61437
SMARCB1 antibody (mAb)	ICC, IF, WB	61649	TEAD1 antibody (pAb)	WB	61643
SMARCC1 / BAF155 antibody (pAb)	ICC, IF, WB	61409	Tet1 antibody (mAb)	ChIP-Seq, ICC, IF, IP, WB	61741
SMARCC2 / BAF170 antibody (pAb)	ICC, IF, WB	61471	Tet1 antibody (pAb)	ChIP, ChIP-Seq, IHC, WB	61443
SMARCE1 antibody (pAb)	WB	61581	AbFlex® TET1 antibody (rAb)	ChIP-Seq	91171
SMC1 phospho Ser957 antibody (pAb)	WB	39631	Tet2 antibody (mAb)	ICC, IF, IP, WB	61389
SMC1-α antibody (pAb)	ChIP-Seq, IP, WB	61067	Tet3 antibody (mAb)	ICC, IF, IP, WB	61743
SMC3 antibody (pAb)	WB	61131	Tet3 antibody (pAb)	ChIP, WB	61395
SMC4 antibody (pAb)	IP, WB	61129	TIEG1 / KLF10 antibody (pAb)	WB	61117
SMRT / NCoR2 antibody (mAb)	WB	61105	TNF antibody (mAb)	ELISA	61601
SMRT / NCoR2 antibody (pAb)	WB	39845	TORC2 antibody (pAb)	IP, WB	39779
Smyd3 antibody (pAb)	IHC, WB	61407	TRIM37 antibody (pAb)	WB	61747
SNAI1 / Snail antibody (mAb)	IHC, WB	61367	TROY/TAJ antibody (pAb)	WB	40937
AbFlex® Snail antibody (rAb)	WB	91255	TWIST antibody (mAb)	ICC, IF, IHC, WB	61097
SND1 antibody (mAb)	ICC, IF, IHC, IP, WB	61473	AbFlex® TWIST antibody (rAb)	ChIP, ChIP-Seq, WB	91323
Sox11 antibody (pAb)	WB	61181	Ubiquitin antibody (mAb)	WB	39741
Sox2 antibody (pAb)	ICC, IF, IHC, WB	39823	AbFlex® Ubiquitin antibody (rAb)	WB	91257
Sox2 antibody (pAb)	ChIP, ChIP-Seq, ICC, IF, WB	39843	Ubn1 antibody (pAb)	WB	39869
SOX6 antibody (pAb)	WB	61749	Uhrf1 antibody (mAb)	ICC, IF, IHC, IP, WB	61341
Sp1 non-phospho antibody (pAb)	WB	39809	Uhrf1 antibody (pAb)	WB	39625
Sp1 phospho Ser101 antibody (pAb)	WB	39757	UTF1 antibody (pAb)	WB	61253
SREBP-1 antibody (mAb)	ChIP, EMSA, IP, WB	39939	WDR5 antibody (pAb)	ChIP, WB	61485
SREBP-2 antibody (mAb)	WB	39941	WRN antibody (mAb)	ICC, IF, IHC, WB	61169
SRF antibody (mAb)	ChIP, ChIP-Seq, IF, IHC, WB	61385	XRCC1 antibody (pAb)	ICC, IF, WB	61371
STAG1 antibody (pAb)	WB	61561	YAP1 antibody (pAb)	IP, WB	61293
STAT1 phospho Ser727 antibody (pAb)	DB, WB	39633	YY1 antibody (pAb)	ChIP-Seq, IHC, IP, WB	61779
STAT2 antibody (pAb)	IP, WB	61651	ZEB1 antibody (pAb)	IP, WB	61119
STAT2 phospho Tyr689 antibody (pAb)	DB, WB	39611			
STAT3 antibody (pAb)	WB	61645			
STAT3 phospho Ser727 antibody (pAb)	DB, WB	39613			
STAT3 phospho Tyr705 antibody (pAb)	DB, WB	39595			
STAT5A antibody (pAb)	IP, WB	61133			
STAT5A/B phospho Ser726/Ser731 antibody (pAb)	DB, WB	39597			
STAT5A/B phospho Tyr694/Tyr699 antibody (pAb)	DB, WB	39617			
Supt16H / Spt16 antibody (pAb)	WB	61441			

The lists above are a small subset of our antibodies, we also offer each in a **convenient sample size**.  
 For an up-to-date list of all available antibodies, please visit [activemotif.com/abs](http://activemotif.com/abs).

## North America

Toll Free: 877 222 9543  
Direct: 760 431 1263  
Fax: 760 431 1351  
E-mail: [tech\\_service@activemotif.com](mailto:tech_service@activemotif.com)

## Europe

UK Free Phone: 0800/169 31 47  
FR Free Phone: 0800/90 99 79  
DE Free Phone: 0800/181 99 10  
Direct: +32 (0)2 653 0001  
Fax: +32 (0)2 653 0050  
E-mail: [eurotech@activemotif.com](mailto:eurotech@activemotif.com)

## Japan

Direct: +81 (0)3 5225 3638  
Fax: +81 (0)3 5261 8733  
E-mail: [japantech@activemotif.com](mailto:japantech@activemotif.com)

## China

Direct: +86 21 20926090  
Cell Phone: 18521362870  
E-mail: [techchina@activemotif.com](mailto:techchina@activemotif.com)

