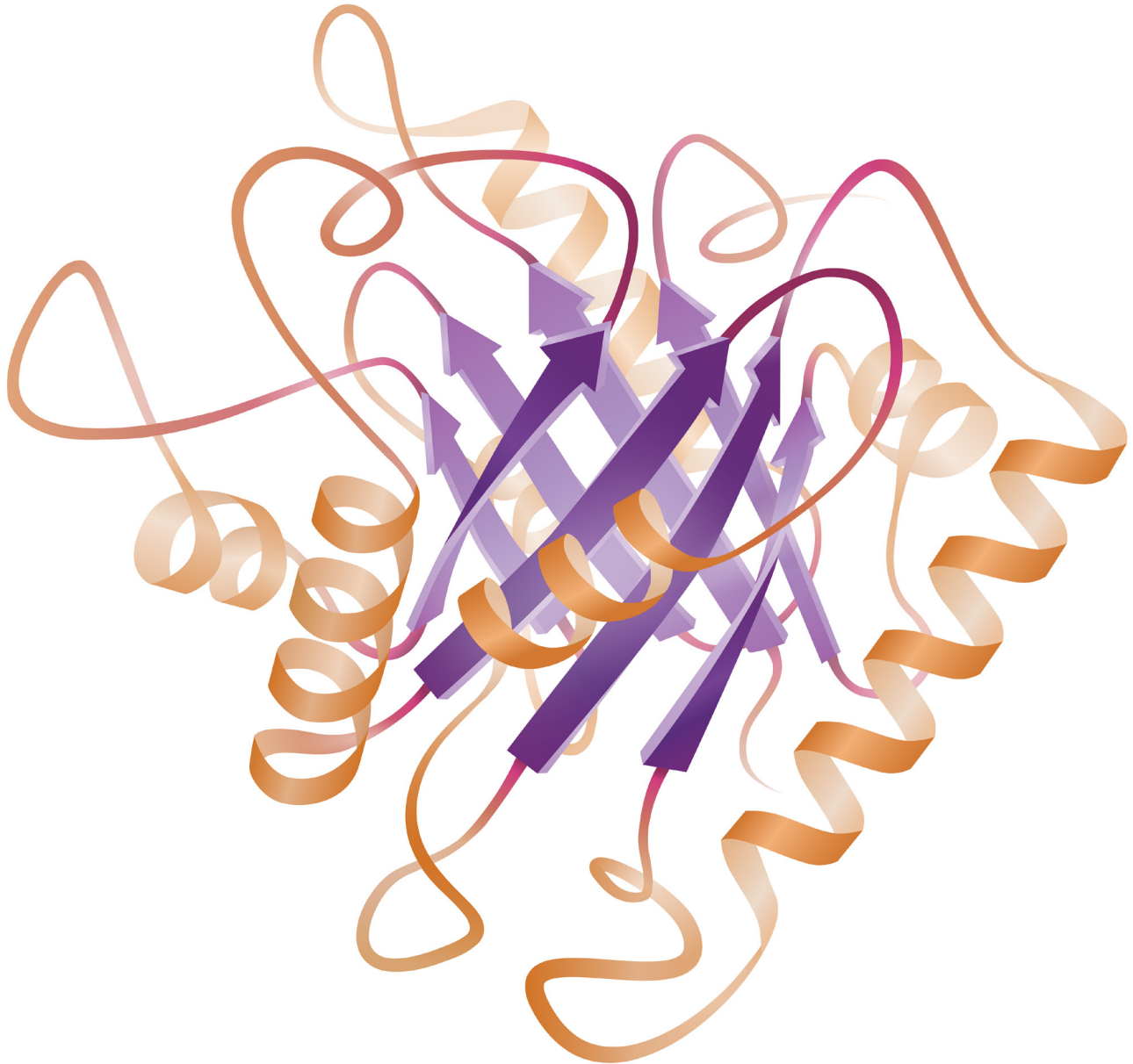


Recombinant Proteins for Epigenetics Research



Choosing the Right Substrate

An important part in the design of enzyme screening assays for epigenetic drug discovery is choosing the correct substrate. For many epigenetic enzymes, screening results are most relevant when the assay most closely simulates physiological conditions. Therefore, choosing biologically relevant substrates such as full-length histones, histone octamers, and nucleosomes will greatly enhance enzyme performance and ultimately, your assay results.

Active Motif offers the most comprehensive collection of recombinant histones, histone octamers, and pre-assembled nucleosomes on the market, giving you the flexibility to choose the best substrate for your assay.

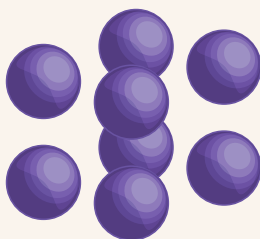
The largest collection of epigenetic substrates

- Unmodified & Modified Histones
- Histone Octamers
- Nucleosomes
- Biotinylated Substrates

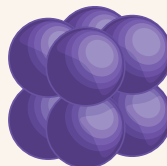
Flexible options for every assay type



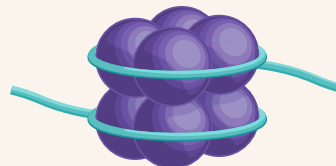
Peptides



Histones



Octamers



Nucleosome



View a complete list at
[activemotif.com/recombhis](https://www.activemotif.com/recombhis)

Recombinant Histones

Active Motif offers a wide variety of full-length recombinant histones that include site- and degree-specific modifications, such as methylation and acetylation. Our modified histones are prepared using patented synthesis technologies to generate histones that most closely mimic native histone substrates. We also offer biotinylated versions of our histone H3 proteins for use in FRET assays and bead-based capture techniques.

Active Motif offers over 150 different recombinant unmodified and modified H1, H2A, H2B, H3 and H4 histones. Use them as stand-alone substrates or assemble to generate complete nucleosomes and oligonucleosomes.

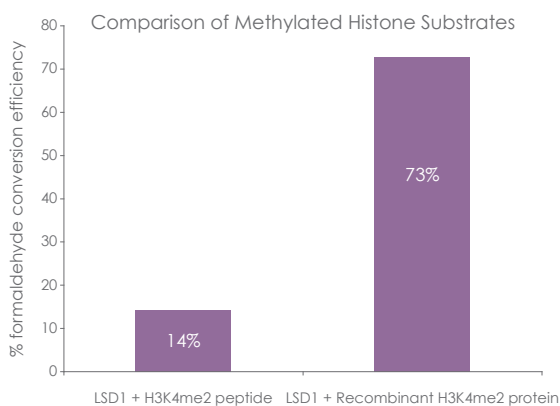


Figure 1: Epigenetic Enzymes are More Active with Full-Length Histone Substrates

Comparison of demethylase efficiency using the Histone Demethylase Assay shows LSD1 demethylates H3K4me2 protein substrate more efficiently than H3K4me2 peptide.

Advantages:

Largest collection of full-length histones

Post-translational modifications (PTMs)

Site- and degree-specific modifications

Oncohistones & histone variants

Biotinylated histones available for FRET

Histone Octamers & Nucleosomes

The structural complexity of chromatin presents a unique challenge to development of enzymatic screening assays for epigenetic drug discovery because of the difficulty in reproducing the chromatin structure *in vitro*. The performance of many epigenetic enzymes in biochemical assays is greatly enhanced when a more physiologically relevant substrate such as an intact nucleosome is available as a substrate.

Active Motif offers a suite of novel, pre-assembled Recombinant Octamers, Mononucleosomes, and Polynucleosomes for use in enzyme activity assays. Additionally, choose from unlabeled or biotin labeled nucleosomes to give you flexibility in your experimental design.

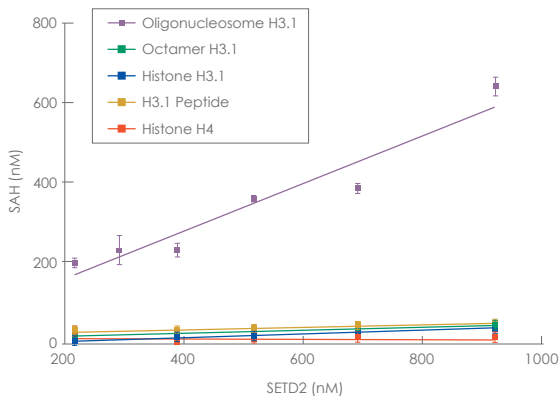


Figure 2: SETD2 Activity Requires Assembled Nucleosomes for its Activity

SETD2 has an absolute requirement for nucleosome substrates.

Comparison of SETD2 activity towards a variety of substrates was measured using an HTRF assay detecting conversion of SAM to SAH. This assay format enables direct comparison of the various substrates but requires more enzyme.

Pre-Assembled Substrates:

Histone Octamers

Mononucleosomes

Oligonucleosomes

Biotinylated versions available

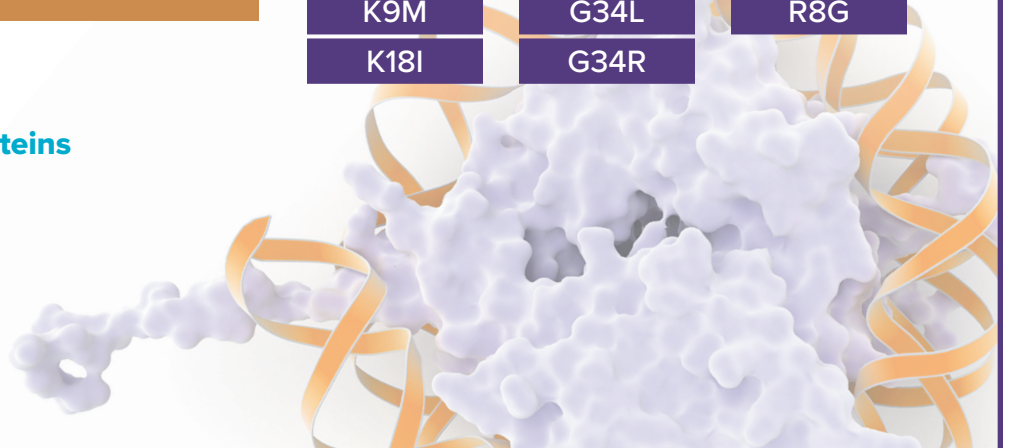
Oncogenic Amino Acid (a.a.) Point Mutations

available as Recombinant Onconucleosomes

K4I	K18M	G34V
K4M	K27M	G34W
K9I	K36M	R8C
K9M	G34L	R8G
K18I	G34R	

View a complete list at

activemotif.com/epiproteins



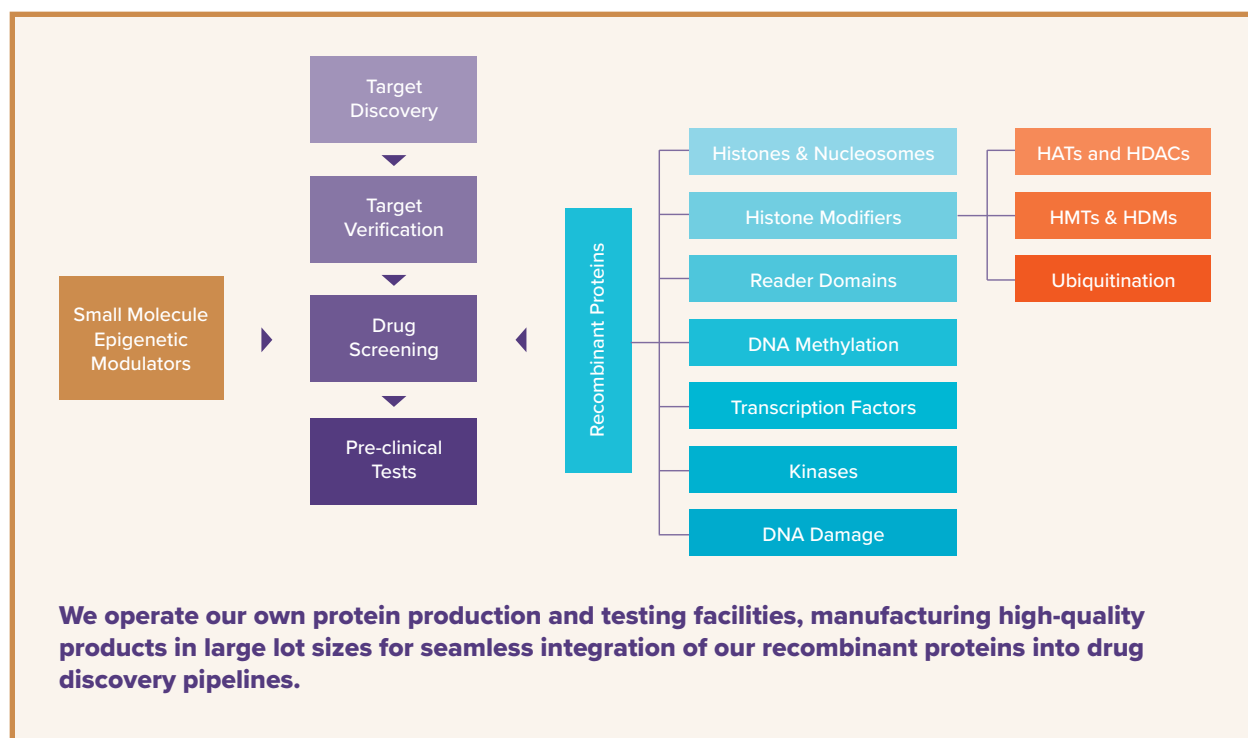
Epigenetic Enzymes & Readers

Active Motif offers a comprehensive portfolio of over 600 ready-to-use purified recombinant epigenetic proteins for use in epigenetic drug discovery research. Our recombinant proteins are manufactured and validated in-house using conventional drug discovery testing methods such as HTRF, mass spectrometry, AlphaLISA, and AlphaScreen.

Our strict quality control measures ensure our proteins meet the highest purity and activity requirements for these types of assays to guarantee peak performance in your research.

Proteins for Epigenetic Drug Discovery Research

- Methyltransferases & Demethylases
- Acetylases & Deacetylases
- Bromodomains
- 1 mg sizes available, ready to ship



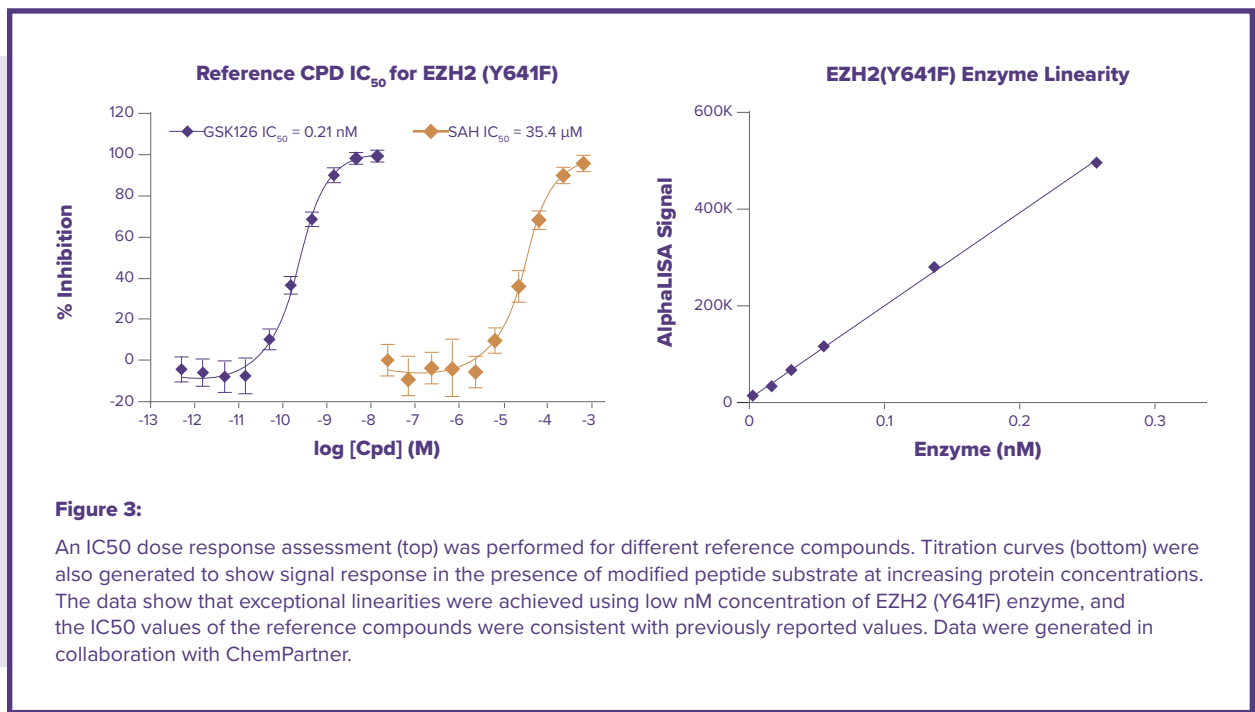
View a complete, up-to-date list of available recombinant proteins at [activemotif.com/proteins](https://www.activemotif.com/proteins)

Methyltransferases & Demethylases

Targeting epigenetic modifiers, in particular methyltransferases and demethylases, has become a primary focus for the development of anti-cancer therapies. In recent years, many novel compounds have been identified that modulate histone and DNA methylation, and several have been developed into therapeutic drugs or moved into clinical trials.

Active Motif offers the largest selection of full-length high quality, robust methyltransferases and demethylases for use in the development of activity assays for drug discovery research. We have produced N-terminal FLAG-tagged HMTs, HDMs, PRMTs, PRDMs, and enzyme complexes for many relevant drug targets, such as NSD2, DOT1L, LSD1, KDM5A, KDM5B, PRC2, and more.

DNA methyltransferases (DNMTs) and oxidation (TET) enzymes are also available for use in studies of DNA methylation.



We provide proteins for the most relevant HMT and HDM drug targets.

Key Products:

KMTs & KDMs

PRMTs & PRDMs

DNMTs

TET Enzymes

Acetyltransferases & Deacetylases

High Quality HAT & HDAC Enzymes

With the FDA approval of histone deacetylase (HDAC) inhibitors as cancer therapies, HDACs and histone acetyltransferases (HATs) have become a focus for drug intervention studies. However, the mechanism of action and specificity of these enzymes continue to be investigated and the pharmacological implications of targeting lysine acetylation as a novel therapeutic strategy is still being explored.

Active Motif provides highly active, high quality HAT & HDAC enzymes to aid in your histone acetylation research. Our HDAC collection includes Class I & II HDACs as well as Class III Sirtuins.

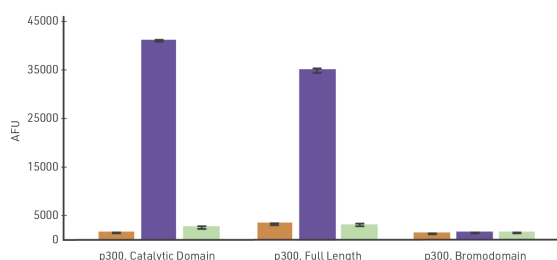


Figure 4:

Activity of p300, Catalytic Domain (Cat. No. 81093), p300, Full Length (Cat. No. 81158), and p300, Bromodomain (Cat. No. 31372) was assessed using the Active Motif HAT Assay Kit (Cat. No. 56100). Equimolar amounts of p300 proteins were incubated with H3 peptide for 3 hours, and developed as per the kit protocol. Anacardic acid, an inhibitor of HAT activity, was used at 15 μ M / well. Graph shows averages \pm SEM of arbitrary fluorescence units (AFU).

Choose From:

p300 full length and catalytic domains

Class I & Class II HDACs

KATs & Sirtuins (SIRTs)

Small Molecule Inhibitors of acetylation & deacetylation



Bulk sizes available. Ready to ship!
[activemotif.com/epiproteins](https://www.activemotif.com/epiproteins)

Reader Domains

Bromodomains for Epigenetic Drug Discovery

Bromodomain (BRD) proteins play an integral role in the regulation of transcription and chromatin remodeling by acting as 'readers' of acetylated histone lysine residues. Because bromodomain proteins have been shown to also regulate transcription of certain oncogenes, they are promising therapeutic targets for cancer.

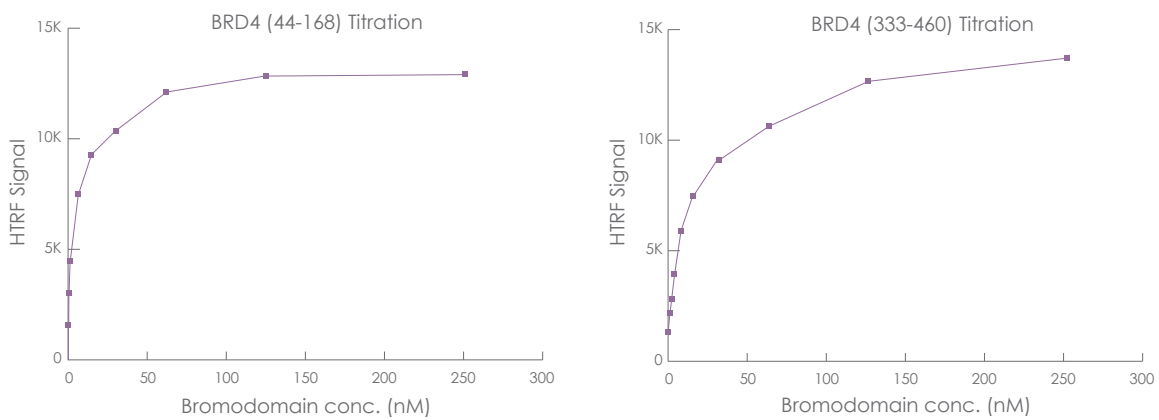


Figure 5: Analysis of activity associated with specific bromodomain regions BRD4 (44-168) (left panel) and BRD4 (333-460) (right panel) bromodomains tested by HTRF.

Assay conditions for bromodomain activity were as follows: 3.3 μ M histone peptide H4K5/8/12/16(tetra-acetyl) was incubated with the protein indicated in reaction buffer containing 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature. Anti-FLAG antibody was used to detect the reaction products.

Active Motif Advantages

- Over 50 reader domains available
- Small molecule inhibitors of Bromodomain containing proteins
- Companion ChIP-Seq validated antibodies for BRD proteins

Additional Drug Discovery Proteins

Proteins for Gene Regulation & Nuclear Biology Research

Active Motif also provides an expansive offering of proteins to study gene regulation and nuclear function including transcription factors, growth factor receptors, nuclear receptors, protein kinases, as well as ubiquitination and neddylation enzymes.

Choose From:

Transcription Factors

Kinases

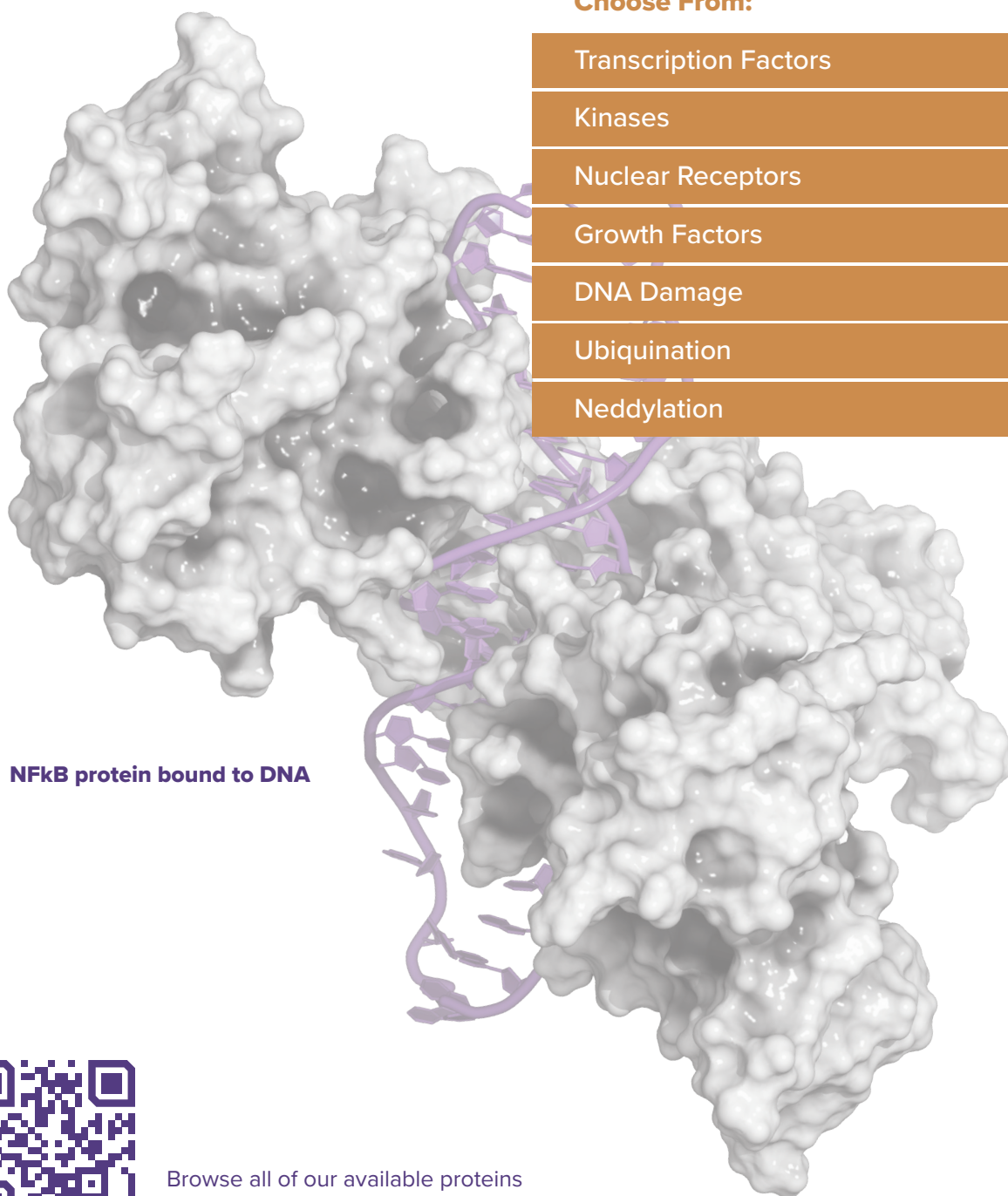
Nuclear Receptors

Growth Factors

DNA Damage

Ubiquitination

Neddylation



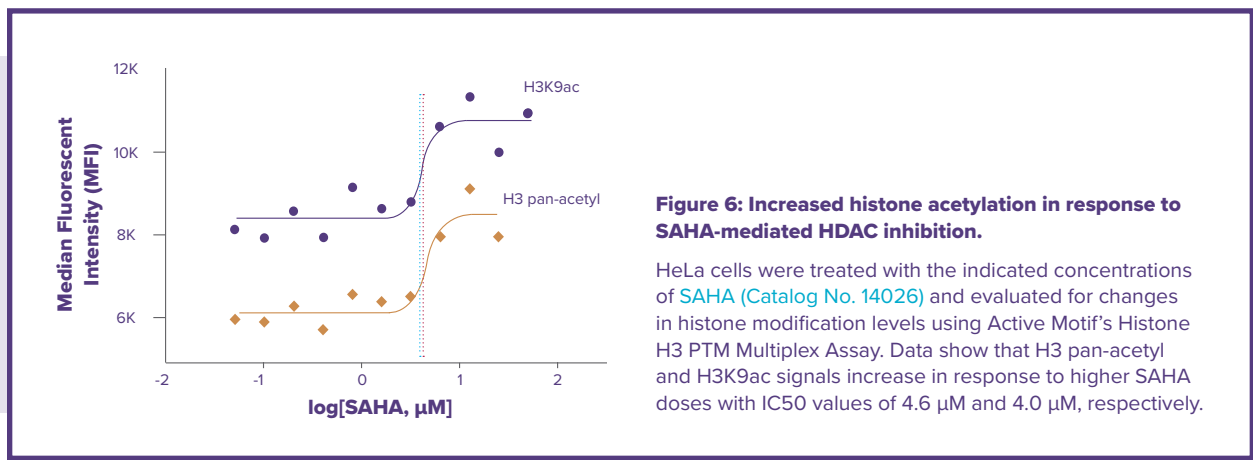
NFkB protein bound to DNA



Browse all of our available proteins
[activemotif.com/epiproteins](https://www.activemotif.com/epiproteins)

Small Molecules to Modulate Activity

Active Motif offers an expanding collection of small molecules (activators and inhibitors) to modulate the activity of proteins that regulate DNA methylation, histones and chromatin. These compounds are ready for use as epigenetic drug discovery tools for lead generation and assay development.



Activators & Inhibitors of:

Lysine Methyltransferases

Lysine Demethylases

Arginine Methyltransferases

Lysine Acetylases

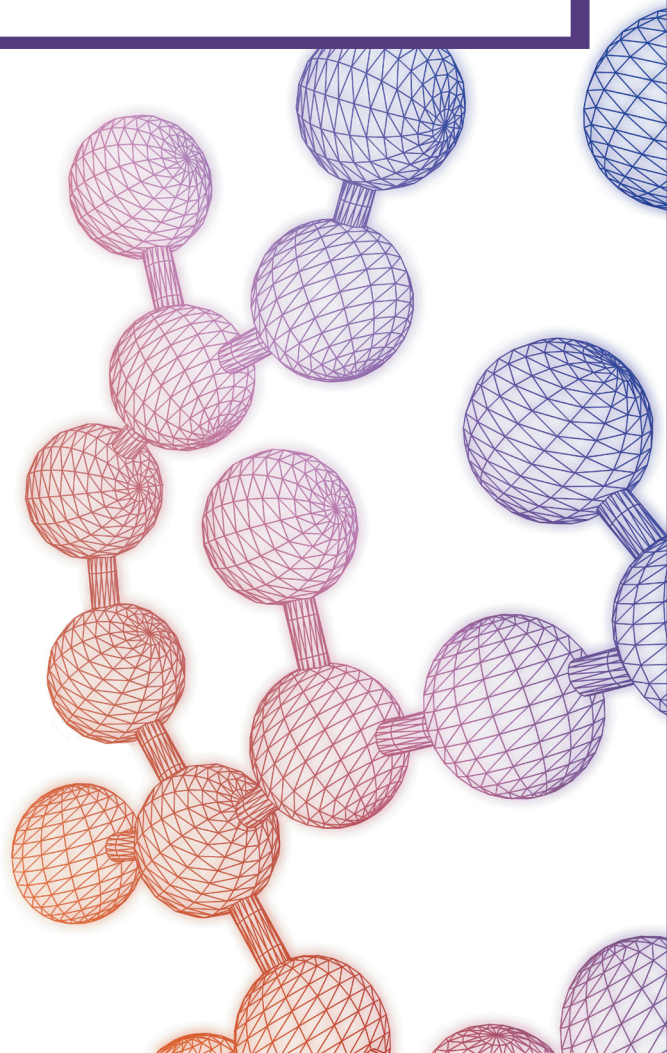
Lysine Deacetylases

Bromodomains

DNA Methylation

For a complete, up-to-date list of available activators and inhibitors, please visit our website at

activemotif.com/smallmol



Ordering Information

Product	Cat. No.	Product	Cat. No.
ABL1 (229-500) protein	81334	BRD2 (71-194) protein, GST-Tag	81149
ABL2 (279-546) protein	81338	BRD2 (BD1+BD2) protein	81435
ACE2 (18-740) protein	81351	BRD3 (24-144) protein	31379
ACF complex	31509	BRD3 (24-144) protein, GST-Tag	81151
AGO1 protein	31522	BRD3 (306-416) protein	31377
AGO2 protein	31486	BRD3 (306-416) protein, GST-Tag	81152
AGO3 protein	31523	BRD3 (BD1+BD2) protein	81436
AKT1 protein	31511	BRD4 (333-460) protein	31446
AKT1 protein	81145	BRD4 (333-460) protein, GST-Tag	81154
AKT2 protein	81146	BRD4 (44-168) protein	31380
AKT3 protein	81147	BRD4 (44-168) protein, GST-Tag	81153
ALDOA protein	81304	BRD4 (44-460) protein	31594
ALDOC protein	81305	BRD4 (44-460) protein, GST-Tag	81155
ALKBH1 protein	81128	BRD7 (129-236) protein	31381
ALKBH1 protein, FLAG-Tag	81308	BRD7 (129-236) protein, GST-Tag	31480
ALKBH2 protein	81129	BRD9 (130-259) protein	31382
ALKBH3 protein	81130	BRD9 (130-259) protein, GST-Tag	31488
ALKBH4 protein	81131	BRDT (21-137) protein	31450
ALKBH5 protein	31589	BRDT (21-137) protein, GST-Tag	81156
ALKBH7 protein	81132	BRDT (257-382), GST-Tag	31483
ALKBH8 protein	81197	BRPF1 (627-746) protein	31375
AMPK Complex (A1+B1+G1) protein	81438	BRPF3 (576-701) protein	31487
AMPK Complex (A1+B1+G2) protein	81437	BTK protein	81083
AMPK Complex (A1+B2+G2) protein	81405	c-Fos protein	31115
AMPK Complex (A2+B2+G1) protein	81406	c-Jun protein	31116
AMPK Complex (A2+B2+G2) protein	81407	c-Myc / MAX complex	81087
APOBEC3A (A3A) protein	81285	c-Myc protein	31117
AS3MT protein	81392	Cas9 protein	81067
ASH1L (2407-2579) protein	31445	CDKN1A protein	81434
AURKA protein	81337	CDKN1B protein	81390
AURKB protein	81352	CDKN2A protein	81404
AXL (470-894) protein	81335	CDKN2C protein	81389
beta-Glucosyltransferase protein	81249	CENPB protein	81201
BirA protein, His-Tag	31517	CHD1 protein	81307
BPTF / FALZ (2791-2911) protein	31447	CHD3 protein	81386
BRCA1 protein	31113	CHD4 protein	81385
BRD1 (556-688) protein	31438	CIAPIN1 protein	81394
BRD2 (344-455) protein	31378	CREBBP (1075-1873) protein	31590
BRD2 (344-455) protein, GST-Tag	81150	CREBBP (1081-1197) protein	31373
BRD2 (71-194) protein	31442	CTBP1 protein	81250

Product	Cat. No.
CTBP2 protein	81251
DAPK3 / ZIPK protein	81432
dCas9 protein, His/AM Tag	81068
dCas9 protein, His/FLAG Tag	81069
DmTet (1634-1986, 2601-2708) protein	81377
DNMT1 protein	31335
DNMT1 protein	31404
DNMT3A (278-432) protein	31541
DNMT3A / DNMT3L complex	31415
DNMT3A protein	31406
DNMT3B (212-358) protein	31542
DNMT3B / DNMT3L complex	31416
DNMT3B protein	31413
DNMT3L protein	31414
DOT1L (1-416) protein	31474
EEF1A1 protein	81219
EEF1A2 protein	81220
EEF1B2 protein	81221
EEF1G protein	81222
EGFR (672-1210) protein	81271
EGFR protein	31165
EGFR protein (672-1210, L858R) protein	81200
EHMT1 (894-1298) protein	31520
EHMT2 (G9a) protein	31410
EHMT2 (G9A)-SET (913-1193) protein	31425
EIF4A1 protein	81216
EIF4A2 protein	81217
EIF4A3 protein	81218
EIF4EBP1 protein	81391
EMG1 protein	81401
Estrogen Receptor α protein	31119
Estrogen Receptor α protein	81276
EZH1 complex	31500
EZH2 protein complex	31337
FAK (409-698) protein	81144
FAK protein	31168
FAK protein	81108
FBPase1 protein	81306
FBXL10 / KDM2B protein	31455
FLT3 (571-993) protein	81381

Product	Cat. No.
FRK (208-505) protein	81357
FTO protein	31572
Fumarase / FH (45-510) protein	81293
Fumarase / FH protein	81292
FXR protein	31120
FYN (2-537) protein	81368
G6PD protein	81299
G6PDH protein	81133
GAMT protein	81393
GLU protein	81298
GNMT protein	81301
GNMT protein	81402
GOT1 (AST1) protein	81297
GR protein	31121
GRK5 protein	81382
GSG2 (470-798) protein	81374
GSK3 β protein	81194
HDAC1 protein	31504
HDAC10 (2-631) protein	81312
HDAC11 (2-347) protein	31538
HDAC2 protein	31505
HDAC3 / NCOR2 complex	31526
HDAC3 / NCOR2 complex, His-Tag	31609
HDAC4 (627-1084) protein	31527
HDAC4 protein	31364
HDAC5 protein	31534
HDAC6 (597-728, deleted mutant) protein	31506
HDAC6 (H230A) protein	31564
HDAC6 protein	31543
HDAC7 (518-991) protein	31535
HDAC8 protein	31536
HDAC8 protein, His-Tag	31566
HDAC9 (604-1066) protein	31537
HER4 / ErbB4 (682-993) protein	81356
HIPK2 (1-640) protein	81355
HIPK3 (163-562) protein	81364
Histone H3K27 Peptide - biotinylated	81048
Histone H3K27ac Peptide - biotinylated	81049
Histone H3K27me1 Peptide - biotinylated	81050
Histone H3K27me2 Peptide - biotinylated	81051

Product	Cat. No.
Histone H3K27me3 Peptide - biotinylated	81052
Histone H3K36 Peptide - biotinylated	81053
Histone H3K36ac Peptide - biotinylated	81054
Histone H3K36me1 Peptide - biotinylated	81055
Histone H3K36me2 Peptide - biotinylated	81056
Histone H3K36me3 Peptide - biotinylated	81057
Histone H3K4 Peptide - biotinylated	81038
Histone H3K4ac Peptide - biotinylated	81039
Histone H3K4me1 Peptide - biotinylated	81040
Histone H3K4me2 Peptide - biotinylated	81041
Histone H3K4me3 Peptide - biotinylated	81042
Histone H3K9 Peptide - biotinylated	81043
Histone H3K9ac Peptide - biotinylated	81044
Histone H3K9me1 Peptide - biotinylated	81045
Histone H3K9me2 Peptide - biotinylated	81046
Histone H3K9me3 Peptide - biotinylated	81047
Histone H4K20 Peptide - biotinylated	81109
Histone H4K20ac Peptide - biotinylated	81110
Histone H4K20me1 Peptide - biotinylated	81111
Histone H4K20me2 Peptide - biotinylated	81112
Histone H4K20me3 Peptide - biotinylated	81113
Histone H1	81126
Histone H1.2	81252
Histone H2A (Human)	31490
Histone H2A.Z/H2B dimer	81168
Histone H2A/H2B dimer	81167
Histone H2B (Human)	31492
Histone H2BFWT	31578
Histone H3 (C110A)	31207
Histone H3.1 (E105K)	81235
Histone H3.1 (E105Q)	81236
Histone H3.1 (E97K)	81234
Histone H3.1 (Human)	31294
Histone H3.1 (K18I)	81239
Histone H3.1 (K18M)	81231
Histone H3.1 (K27M)	81233
Histone H3.1 (K4I)	81228
Histone H3.1 (K4M)	81226
Histone H3.1 (K9I)	81229
Histone H3.1 (K9M)	81227

Product	Cat. No.
Histone H3.1 (R26C)	81232
Histone H3.1 (R26H)	81241
Histone H3.1 (R26P)	81240
Histone H3.1 (R8G)	81230
Histone H3.1 biotinylated (Human)	31296
Histone H3.1t	81238
Histone H3.2	81237
Histone H3.2 biotinylated (Human)	31271
Histone H3.3 (G34R)	31549
Histone H3.3 (G34V)	31550
Histone H3.3 (G34W)	31551
Histone H3.3 (Human)	31295
Histone H3.3 (K18I)	81248
Histone H3.3 (K18M)	81247
Histone H3.3 (K27M)	31552
Histone H3.3 (K36M)	31553
Histone H3.3 (K4I)	81244
Histone H3.3 (K4M)	81243
Histone H3.3 (K9I)	81246
Histone H3.3 (K9M)	81245
Histone H3.3 (R8C)	81242
Histone H3.3 biotinylated (Human)	31297
Histone H3/H4 tetramer	81169
Histone H3K14ac (EPL)	31254
Histone H3K14me1 (MLA)	31256
Histone H3K14me2 (MLA)	31257
Histone H3K14me3 (MLA)	31258
Histone H3K18ac (EPL)	31273
Histone H3K18me1 (MLA)	31259
Histone H3K18me2 (MLA)	31260
Histone H3K18me3 (MLA)	31261
Histone H3K23ac (EPL)	31255
Histone H3K23me1 (MLA)	31262
Histone H3K23me2 (MLA)	31263
Histone H3K23me3 (MLA)	31264
Histone H3K27me1 (MLA)	31621
Histone H3K27me2 (MLA)	31597
Histone H3K27me3 (MLA)	31579
Histone H3K36me1 (MLA)	31622
Histone H3K36me2 (MLA)	31598

Product	Cat. No.
Histone H3K36me3 (MLA)	31580
Histone H3K4ac (EPL)	31275
Histone H3K4me1 (EPL)	31287
Histone H3K4me1 (MLA)	31619
Histone H3K4me1 biotinylated (EPL)	31284
Histone H3K4me2 (EPL)	31277
Histone H3K4me2 (MLA)	31595
Histone H3K4me2 biotinylated (EPL)	31283
Histone H3K4me3 (EPL)	31278
Histone H3K4me3 (MLA)	31600
Histone H3K4me3 biotinylated (EPL)	31282
Histone H3K79me1 (MLA)	31220
Histone H3K79me2 (MLA)	31599
Histone H3K79me3 (MLA)	31581
Histone H3K9ac (EPL)	31253
Histone H3K9me1 (EPL)	31281
Histone H3K9me1 (MLA)	31620
Histone H3K9me1 biotinylated (EPL)	31286
Histone H3K9me2 (EPL)	31280
Histone H3K9me2 (MLA)	31596
Histone H3K9me3 (EPL)	31279
Histone H3K9me3 (MLA)	31601
Histone H3K9me3 biotinylated (EPL)	31285
Histone H3R8me2a (asymmetric) (EPL)	31276
Histone H3S10ph (EPL)	31272
Histone H3T3ph (EPL)	31274
Histone H4, His-Tag (Human)	31493
Histone H4K16me1 (MLA)	31268
Histone H4K16me2 (MLA)	31269
Histone H4K16me3 (MLA)	31270
Histone H4K20me3 (MLA)	31226
Histone H4K5me1 (MLA)	31265
Histone H4K5me2 (MLA)	31266
Histone H4K5me3 (MLA)	31267
Histone Octamer (H3.1)	31470
Histone Octamer (H3.1) - biotinylated	31471
Histone Octamer (H3.3)	31472
Histone Octamer (H3.3) - biotinylated	31473
Histone TH2B	31577
hnRNPA2B1 protein	31607

Product	Cat. No.
IDE protein	81303
IDH1 (R132C) protein	31613
IDH1 (R132H) protein	31614
IDH1 protein	31610
IDH2 (R140K) protein	31616
IDH2 (R140Q) protein	31617
IDH2 (R172K) protein	31618
IDH2 protein	31615
IDO1 protein	81031
IDO2 (15-420) protein	31587
IGF1R (763-931) protein	81358
IGF1R (960-1367) protein	81360
IGF2BP1 protein	81289
IGF2BP2 protein	81290
IGF2BP3 protein	81291
IKK β protein	81066
IKK ϵ protein	81117
INSR (1011-1382) protein	81361
INSR (999-1362) protein	81359
IRF3 protein	31544
ITK (352-620) protein	81362
JAK1 (438-1154) protein	81369
JAK1 (866-1154) protein	81370
JAK2 (532-1132, V617F) protein	81283
JAK2 (532-1132) protein	81311
JAK3 (781-1124) protein	81373
JARID1A / KDM5A protein	31431
JARID1B / KDM5B (2-751) protein	31518
JARID1B / KDM5B protein	31432
JARID1C / KDM5C protein	31433
JHDM1D-I protein	31464
JMJD1A / KDM3A protein	31456
JMJD1B / KDM3B protein	31429
JMJD2A / KDM4A protein	31457
JMJD2B / KDM4B protein	31501
JMJD2C / KDM4C protein	31458
JMJD2D / KDM4D protein	31459
JMJD3 / KDM6B (1043-1682) protein	31519
KAT1 / HAT1 protein	81274
KAT2A (GCN5) protein	31591

Product	Cat. No.
KAT2B / PCAF (715-829) protein	31370
KAT2B / PCAF protein	81142
KAT5 protein	81275
KAT6A / MOZ (488-778) protein	81223
KAT6B / MORF (718-1008) protein	81224
KAT7 protein	31489
KAT8 / MYST1 protein	81225
KDM1B / LSD2 protein	31479
KDM2A / FBXL11 protein	31485
KDM6C protein	31575
KDM7A protein	31576
KDR / VEGFR2 (789-1356) protein	81353
Ketohexokinase protein	81300
KIT / CD117 (546-976) protein	81371
KMT2A (MLL1) complex	31423
KMT2A (MLL1)-SET protein	31419
KMT2B (MLL4) complex	31499
KMT2B (MLL4)-SET protein	31422
KMT2C (MLL3) complex	31478
KMT2D (MLL2) complex	31498
KMT2D (MLL2)-SET protein	31420
LATS1 protein	81209
LCK protein	81384
LSD1 / KDM1A protein	31426
LXR α protein	31122
LXR β protein	31123
LYN protein	81367
MAOA protein	31502
MAOB protein	31503
MAP2K1 (MEK1) protein	81199
MAP2K2 protein	81332
MAP3K5 (660-978) protein	81333
MAP3K8 (30-397) / COT protein	81380
MAPK3 (ERK1) protein	81157
MAPKAPK3 protein	81366
MAX protein	81017
MAX protein, His-Tag	81026
MDH1 protein	81295
MDH2 (25-338) protein	81296
MER (528-999) protein	81375

Product	Cat. No.
METTL1 / WDR4 complex	81206
METTL1 protein	81022
METTL1 protein, His-Tag	81058
METTL10 protein	81212
METTL13 protein	81024
METTL14 protein	31568
METTL15 protein	81084
METTL16 protein	81085
METTL18 protein	81030
METTL19 (TRMT44) protein	81088
METTL21B protein	81207
METTL21D (VCPKMT) protein	81210
METTL22 protein	81092
METTL25 protein	81089
METTL2A protein	81027
METTL3 / METTL14 complex	31570
METTL3 protein	31567
METTL4 protein	81177
METTL5 protein	81211
METTL6 protein	81023
METTL6 protein, His-Tag	81060
METTL7A protein	81178
METTL7B protein	81208
METTL8 protein	81028
METTL8 protein, GST-Tag	81106
MINK1 (1-320) protein	81417
MLLT1 / ENL (1-148) protein	81098
MLLT3 / AF9 (1-138) protein	81124
MLLT3 / AF9 (1-138) protein, His/FLAG Tag	81192
Mononucleosomes (H1.2)	81272
Mononucleosomes (H1.2) - biotin	81273
Mononucleosomes (H2A.Bbd) - biotinylated	31556
Mononucleosomes (H2A.X)	81125
Mononucleosomes (H2A.X) - biotinylated	31582
Mononucleosomes (H2A.Z)	81072
Mononucleosomes (H2A.Z) - biotinylated	31583
Mononucleosomes (H3.1)	81070
Mononucleosomes (H3.1) - biotinylated	31467
Mononucleosomes (H3.3)	81071
Mononucleosomes (H3.3) - biotinylated	31469

Product	Cat. No.
Mononucleosomes (TH2B) - biotinylated	31557
Mononucleosomes H3.1 (K18I)	81260
Mononucleosomes H3.1 (K18I) - biotin	81268
Mononucleosomes H3.1 (K18M)	81255
Mononucleosomes H3.1 (K18M) - biotin	81263
Mononucleosomes H3.1 (K27M)	81256
Mononucleosomes H3.1 (K27M) - biotin	81264
Mononucleosomes H3.1 (K4I)	81257
Mononucleosomes H3.1 (K4I) - biotin	81265
Mononucleosomes H3.1 (K4M)	81253
Mononucleosomes H3.1 (K4M) - biotin	81261
Mononucleosomes H3.1 (K9I)	81258
Mononucleosomes H3.1 (K9I) - biotin	81266
Mononucleosomes H3.1 (K9M)	81254
Mononucleosomes H3.1 (K9M) - biotin	81262
Mononucleosomes H3.1 (R26C) - biotin	81324
Mononucleosomes H3.1 (R26P) - biotin	81325
Mononucleosomes H3.1 (R8G)	81259
Mononucleosomes H3.1 (R8G) - biotin	81267
Mononucleosomes H3.2	81345
Mononucleosomes H3.2 - biotin	81326
Mononucleosomes H3.3 (G34R) - biotin	81009
Mononucleosomes H3.3 (G34V) - biotin	81010
Mononucleosomes H3.3 (K18I)	81346
Mononucleosomes H3.3 (K18I) - biotin	81327
Mononucleosomes H3.3 (K18M)	81270
Mononucleosomes H3.3 (K18M) - biotin	81288
Mononucleosomes H3.3 (K27M) - biotin	81006
Mononucleosomes H3.3 (K4I)	81348
Mononucleosomes H3.3 (K4I) - biotin	81328
Mononucleosomes H3.3 (K4M)	81347
Mononucleosomes H3.3 (K9I)	81350
Mononucleosomes H3.3 (K9I) - biotin	81330
Mononucleosomes H3.3 (K9M)	81349
Mononucleosomes H3.3 (K9M) - biotin	81329
Mononucleosomes H3.3 (R8C)	81269
Mononucleosomes H3.3 (R8C) - biotin	81287
Mononucleosomes H3K14ac (EPL) - biotin	81001
Mononucleosomes H3K27ac	81077
Mononucleosomes H3K27ac - biotin	81002

Product	Cat. No.
Mononucleosomes H3K27me3 (MLA)	81134
Mononucleosomes H3K27me3 (MLA) - biotin	81135
Mononucleosomes H3K36me3 (MLA)	81116
Mononucleosomes H3K36me3 (MLA) - biotin	81076
Mononucleosomes H3K4me1 (EPL) - biotinylated	31585
Mononucleosomes H3K4me2 (EPL)	81073
Mononucleosomes H3K4me2 (EPL) - biotin	81074
Mononucleosomes H3K4me3 (EPL) - biotinylated	31584
Mononucleosomes H3K4me3/H3K27ac - biotin	81004
Mononucleosomes H3K9ac (EPL)	81075
Mononucleosomes H3K9ac (EPL) - biotin	81018
Mononucleosomes H3K9me3 (EPL)	31586
Mononucleosomes H3K9me3 (EPL) - biotinylated	31555
Mononucleosomes H3R2/8/17 citrul (EPL)	81165
Mononucleosomes H3S10ph (EPL)	81163
Mononucleosomes H3S10ph (EPL) - biotin	81164
MRCKa / CDC42BPA (1-473) protein	81396
MRCKb / CDC42BPB (1-473) protein	81397
MST1 protein	31355
MST2 protein	81414
Myelin Basic Protein, dephosphorylated	31314
NAP1L1 protein	31508
NAT10 protein	81376
NEK2 protein	81413
NEK7 protein	81410
NFκB p50 protein	81310
NFκB p65 protein	31302
NFκB1 p105 protein	81143
NFκB1 p50 (1-434) protein	81032
NFκB3 (RELA / p65) protein	81086
NgTet1 (1-321) protein	81148
NLK protein	81415
NONO protein	31539
NSD1-SET protein	31475
NSD2 (E1099K) protein	31546
NSD2 (MMSET) protein	31453
NSD2 (MMSET)-SET protein	31476
NSD2-SET (E1099K) protein	31574
NSD3 (WHSC1L1)-SET protein	31477
NSUN1 protein	81170

Product	Cat. No.
NSUN2 protein	81171
NSUN3 protein	81183
NSUN5 protein	81173
NSUN6 protein	81174
NTRK1 (440-796) protein	81411
NTRK3 (454-825) protein	81383
NUAK1 (ARK5) protein	81331
OGT protein	31524
p300 (1041-1161) protein	31372
p300 protein	81158
p300 protein	81858
p300 protein, catalytic domain	81093
p53 (TP53) protein	81091
p53 protein	31103
p53 protein	31465
PAK4 protein	81416
PARP1 protein	81037
PBRM1 (613-734) protein	31385
PD-L1 / CD274 (19-239) protein	81387
PDCD1 / PD1 (25-167) protein	81388
PDGFRA (550-1089) protein	81430
PDGFRB (557-1106) protein	81420
PDK1 protein, GST-Tag	81213
PEL1 protein	81313
PHD1 (EGLN2) protein	81064
PHD2 (EGLN1) protein	81065
PHD3 (EGLN3) protein	81033
PHD3 (EGLN3) protein, FLAG-Tag	81215
PHF8 protein	31435
PHKG2 protein	81418
PIK3R1 protein	81097
PIM2 protein	81365
PKM2 protein	81302
PKN2 protein	81431
PLK1 protein	81429
PLK3 (57-340) protein	81372
PLZF protein	31545
Polynucleosomes (H3.1)	31466
Polynucleosomes (H3.3)	31468
Polynucleosomes H3.3 (G34R)	31559

Product	Cat. No.
Polynucleosomes H3.3 (G34V)	31560
Polynucleosomes H3.3 (G34W)	31561
Polynucleosomes H3.3 (K36M)	31563
PPAR α protein	31125
PPAR β (δ) protein	31126
PPAR γ protein	31127
PPAR γ protein	81214
PRC2 complex	31387
PRC2 EZH2 (A677G) complex	31391
PRC2 EZH2 (Y641C) complex	31389
PRC2 EZH2 (Y641F) complex	31388
PRC2 EZH2 (Y641N) complex	31390
PRDM1 protein	81277
PRDM10 protein	31396
PRDM11 protein	81278
PRDM14 protein	31397
PRDM4 protein	81395
PRDM5 protein	31494
PRDM6 protein	31495
PRDM9 (191-414) protein	31510
PRDM9 (191-415) protein	81118
PRKAA1 complex	81340
PRKCD protein	81422
PRKCE protein	81421
PRKCH protein	81426
PRKCI protein	81424
PRKCZ protein	81425
PRKD2 protein	81419
PRKG1 protein	81423
PRMT1 protein	31411
PRMT2 protein	31392
PRMT3 protein	31412
PRMT4 (CARM1) protein	81107
PRMT5 / MEP50 complex	31521
PRMT5 protein	31393
PRMT6 protein	31394
PRMT7 protein	31395
PTEN protein	81403
PTK6 protein	81339
PTPN1 (1-321) protein	81035

Product	Cat. No.
PTPN1 protein	81034
PTPN2 protein	31592
PXR protein	31144
RAR- α protein	31130
RNA Pol II - CTD protein	81036
RPS6KA1/RSK1 protein	81398
RPS6KA2/RSK3 protein	81408
RPS6KA3/MAPKAPK1B protein	81363
RPS6KA4/RSK4 protein	81399
RPS6KB1 protein	81409
RSBN1 / KDM9 protein	81309
RXR-LBD protein	31135
RXR-LBD protein	31365
RXR- α (RXRA) protein	81082
RXR- β protein	31134
SARS-CoV-2 3C-Like Proteinase (NSP5), GST-Tag	81321
SARS-CoV-2 3C-Like Proteinase (NSP5), His-Tag	81320
SARS-CoV-2 NSP1 protein	81314
SARS-CoV-2 NSP10 / NSP16 complex	81319
SARS-CoV-2 NSP10 protein	81317
SARS-CoV-2 NSP12 / RdRp protein	81354
SARS-CoV-2 NSP16 protein	81318
SARS-CoV-2 NSP2 protein	81323
SARS-CoV-2 NSP7 protein	81315
SARS-CoV-2 NSP8 protein	81316
SARS-CoV-2 NSP9 protein	81322
SARS-CoV-2 Spike protein, S1 (RBD aa319-541)	81343
SARS-CoV-2 Spike protein, S1 (RBD aa319-589)	81344
SETD1A-SET complex	81341
SETD1B-SET complex	81342
SETD2 (1392-2564) protein	31399
SETD2 (1418-1714) protein	31358
SETD3 protein	81279
SETD4 protein	81280
SETD5 (185-416) protein	81281
SETD6 protein	81282
SETD7 protein	31496
SETD8 protein	31427
SETDB1 protein	31452
SETMAR protein	31454

Product	Cat. No.
SIRT1 (193-741) protein	31533
SIRT2 (50-356) protein	31528
SIRT3 (102-399) protein	31529
SIRT4 (25-314) protein	31530
SIRT5 protein	31531
SIRT6 protein	31532
SMARCA2 (636-1131) protein	81439
SMARCA2 / BRM (1367-1511) protein	31449
SMARCA2 / BRM (1367-1511), GST-Tag	31481
SMARCA4 (658-1328) protein	81440
SMARCA4 / BRG1 (1448-1569) protein	31401
SMARCA4 / BRG1 (1448-1569), GST-Tag	31482
SMYD1 protein	31405
SMYD2 protein	31497
SMYD3 protein	31407
SMYD4 protein	31408
SMYD5 protein	31409
SORD protein	81294
Sortase (2A.9) protein	13112
Sortase A5 protein	13101
SP1 protein	81181
SRC protein	81115
SSRP1 / FACT p80 protein	81094
STAT3 protein	81095
STK10 / LOK (1-348) protein	81378
STK11 / LKB1 protein	81379
SUV39H1 (82-412) protein	81020
SUV39H2 protein	81012
SUV420H1 (2-387) protein	81021
SUV420H2 (2-281) protein	81014
SUV420H2 protein	81013
TAF1 (1522-1656) protein	31439
TBP protein	81114
Tet1 (1418-2136) protein	31417
TET2 (1129-2002) protein	31418
Tet3 (824-1795) protein	31421
THR β protein	81120
TMEM173 (STING) (149-379) protein	81182
TRIM24 (862-980) protein	31368
TRIM28 (624-811) protein	31441

Product	Cat. No.
TRIM33 (959-1069) protein	31367
TRIT1 (48-467) protein	81190
TRMT61A protein	81433
TRβ1 protein	31139
TSSK2 protein	81412
UHRF1 protein	81015
USP7 protein	31525
UTX / KDM6A protein	31460
VRK1 protein	81096
vSET (A612L) protein	31402
WNK2 (1-489) protein	81428
WTAP protein	31571
YEATS2 (200-340) protein	81099
YEATS4 (GAS41) protein	81193
YES1 protein	81427
YTHDC1 (325-502) protein	81100
YTHDC1 protein	81176
YTHDC2 (1279-1429) protein	81101
YTHDC2 protein	81198
YTHDF1 (380-533) protein	81102
YTHDF1 protein	31608
YTHDF2 (401-554) protein	81103
YTHDF2 protein	31573
YTHDF3 (407-560) protein	81104
YTHDF3 protein	81204
YY1 protein	81119



Take advantage
of Active Motif
Epigenetics
Resources!



Epigenetics 101: Complete Guide to Understanding Epigenetics

Take a deep dive into epigenetics in our comprehensive learning center.



Epigenetics Blog

Follow our posts about the latest trends in epigenetics research.



eBooks

Stay up to date with the latest eBooks from our epigenetics experts.



Scientific Posters

Download from a collection of our scientific posters.



Product Literature

Download educational guides, product brochures, and references lists.



Publications

Search publications that have used our products and services.



Webinars

Watch the latest tips and techniques from our epigenetics experts.



Videos

Browse our video collection and become an epigenetics expert.



Chromatin Clubs & Epigenetics Events

Keep connected to the epigenetics research community with chromatin clubs, online seminars, virtual conferences, and more epigenetics events.

Check out the chromatin club events that we have scheduled or contact us to request support for a chromatin club you are planning.

North America

Toll Free: 877 222 9543
Direct: 760 431 1263
Fax: 760 431 1351
E-mail: tech_service@activemotif.com

Japan

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

Europe

Direct: +32 (0)2 653 0001
Fax: +32 (0)2 653 0050
E-mail: eurotech@activemotif.com

China

Direct: +86 21 20926090
Cell Phone: 18521362870
E-mail: techchina@activemotif.com

