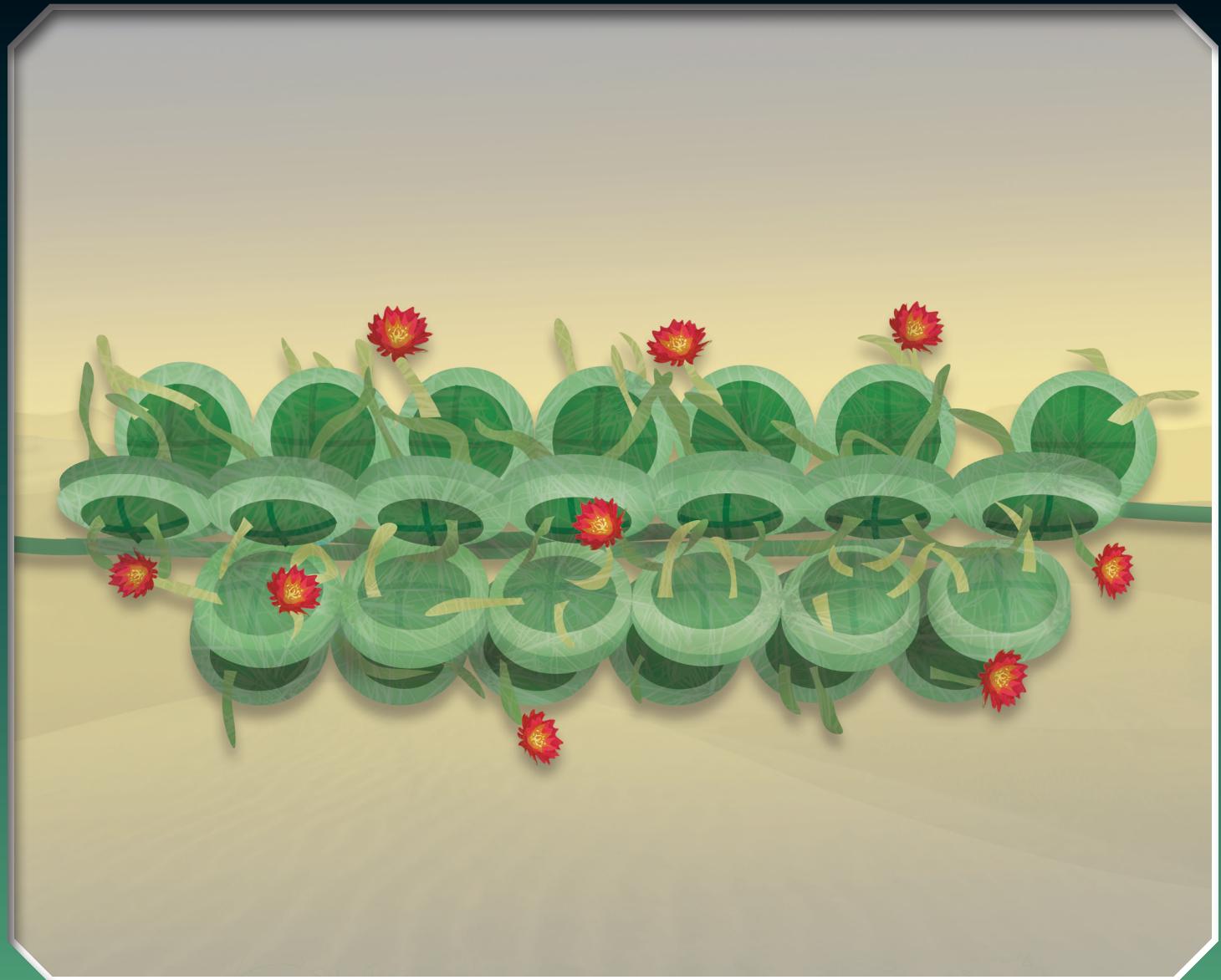




# EPIGENETICS

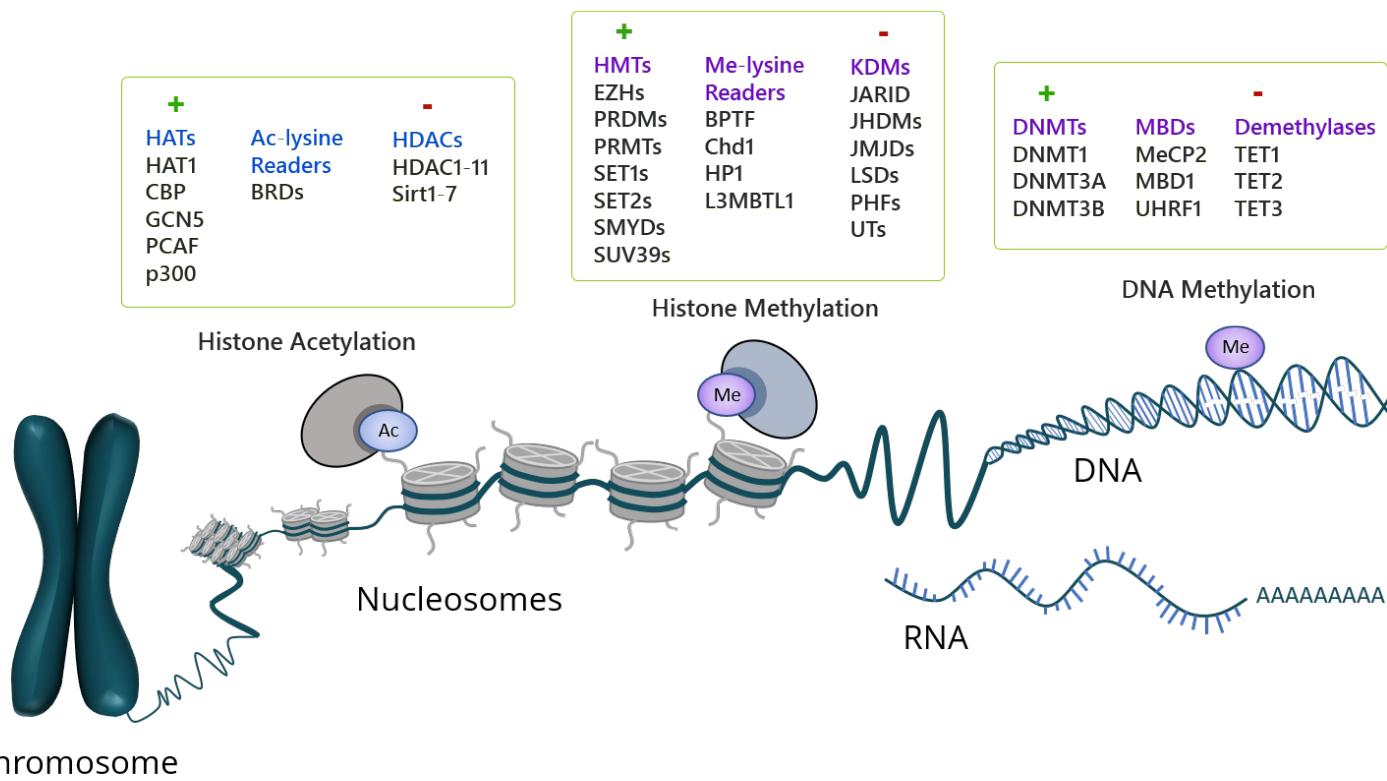
## DRUG DISCOVERY SOLUTIONS

Bioactive Recombinant Enzymes | Assay Kits | Screening & Profiling | Custom Services



# Gene Control Without Gene Editing

The field of epigenetics refers to the control or modulation of gene expression that produces a specific phenotype or trait without directly changing DNA sequences. Epigenetic control can occur by modifying DNA, RNA, or by modulating the proteins that bind to DNA, particularly histone proteins. Epigenetic modifications start early in embryonic development and continue as organisms age and are strongly influenced by environmental factors. Epigenetic mechanisms are now important therapeutic drug targets spanning several disease areas, including multiple cancers, myelodysplastic syndromes, autoimmunity, neurodegeneration, and more.



## Chromosome

Epigenetic control of DNA expression can occur through several mechanisms. Some primary mechanisms include DNA methylation, histone acetylation and histone methylation. The process of addition and removal of these functional groups involves an array of proteins, many of which are indicated in the figure above. These modifications alter chromatin structure and DNA accessibility, thus either preventing or allowing promoter and enhancer regions to be recognized by transcription factors to initiate gene transcription. In addition, there are families of proteins that recognize and bind to these modifications, known as acetyl-lysine readers or methyl-lysine readers, further exerting epigenetic control. Histone can also be further modified by a large number of post-translational modifications, including phosphorylation, ubiquitination, sumoylation, ADP-ribosylation, citrullination, O-GlcNAcylation, and others. In addition to DNA and histones, RNA modifications can also influence translation, resulting in distinct phenotypes, including diseases that are associated with specific RNA modifications.

Our products and services include novel assay kits and bioactive recombinant proteins optimized for drug discovery. The portfolio includes complete sets of specific protein families not found elsewhere. Read on to learn how our solutions can solve your challenges in epigenetic drug discovery.

# Our Advantages



## Produced In-House

- Made in the USA at our San Diego, CA laboratory
- Get customized, personal support directly from the source



## Committed to Excellence

- ISO 9001:2015-certified Quality Management System
- Lot-specific quality testing



## Protein Expression & Screening Services

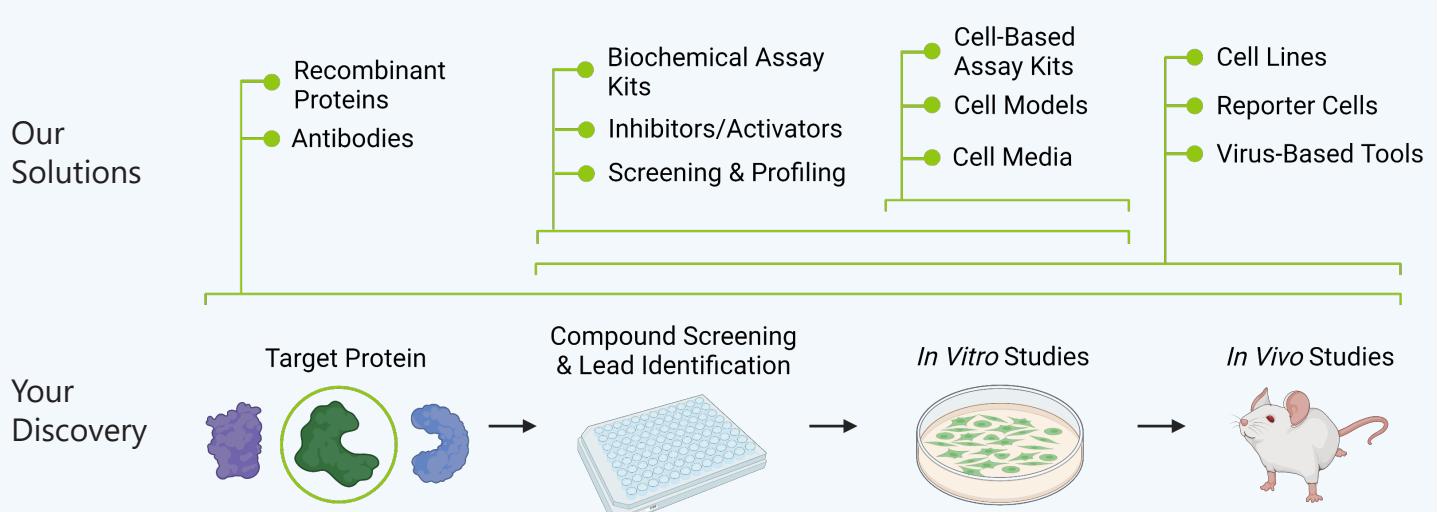
- Screening & Profiling: 110+ biochemical assays to screen your compounds of interest
- Custom protein expression: choice of tags, labeling, expression system, and QC



## Expansive Portfolio

- Assay-ready recombinant proteins: 250+ active epigenetic proteins including complexes
- 175+ inhibitor screening assays
- Substrates - histones, octomers, & nucleosomes, including biotinylated versions
- Full HDAC and sirtuin suite
- Most extensive DNMT offering on the market

## How we drive drug discovery



Created with BioRender.com

# Bioactive Recombinant Proteins

BPS Bioscience specializes in manufacturing bioactive enzymes and protein complexes, offering over 2000 protein products critical for drug discovery. Our epigenetics portfolio consists of over 250 proteins covering complete families of enzymes involved in acetylation and deacetylation, methylation and demethylation, and RNA modification, as well as Bromodomain proteins (acetyl-lysine readers) and methyl-lysine readers.

Protein and peptide substrates, ligands, and small molecule inhibitors are also available.

Protein options include:



## HiP™ Proteins (High Purity, Low Aggregation)

- >90% purity
- Verified via HPLC to guarantee <10% aggregation
- Ideal for studying enzyme activity, ligand binding, protein folding, and more

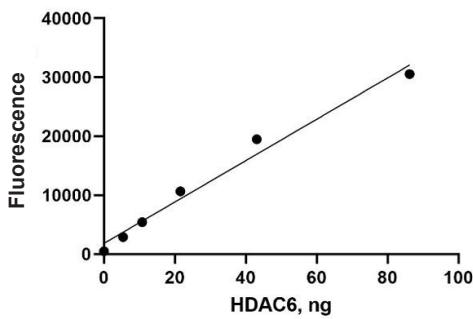
## Bulk Production and Customization

Product sizing, formulations, tags, mutations, and activity testing can be customized to meet your research needs.

## Featuring

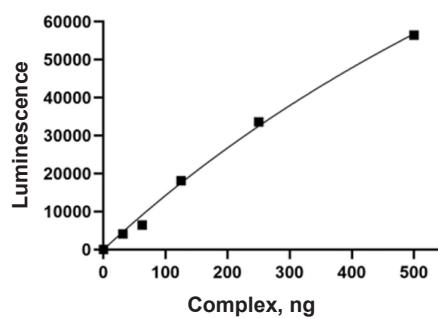
- Active complexes, including DNMT3A/DNMT3L, EZH2/EED/SUZ12/RbAp48/AEBP2, METTL3/METTL14, and more
- Recombinant and native Histone proteins, including tetramers and octamers, with or without tags
- Small molecule inhibitors to establish positive controls for enzyme inhibition
- Mutant and wild-type versions for EZH2, Histone H3, and HDAC6.

### HDAC6 Specific Activity



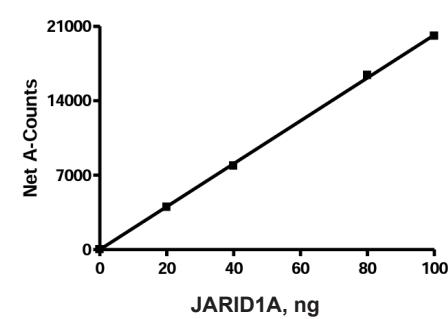
Measurement of HDAC6 activity using a fluorogenic substrate, according to BPS Bioscience #50076, HDAC6 Fluorogenic Assay Kit.

### EZH2/EED/SUZ12/RbAp48/AEBP2 Specific Activity



Measurement of EZH2/EED/SUZ12/RbAp48/AEBP2 complex activity according to BPS Bioscience #52009L, EZH2 Chemiluminescent Assay Kit.

### JARID1A Specific Activity



Measurement of JARID1A demethylation activity using an AlphaLISA® -based assay according to BPS Bioscience #50510, JARID1A Homogeneous Assay Kit.

# Biochemical Assay Kits

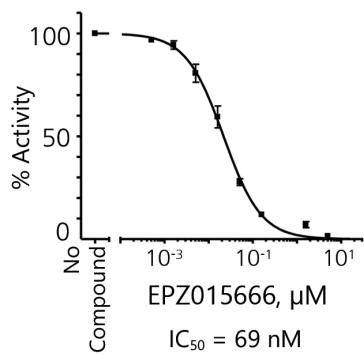
Our 175+ epigenetic assay kits enable your compound and biotherapeutic drug discovery by providing specific, highly sensitive analysis in a variety of formats. Many of our kits are provided as homogenous assays, with no wash steps, for efficient high throughput screening.

Compare the benefits of our assay formats:

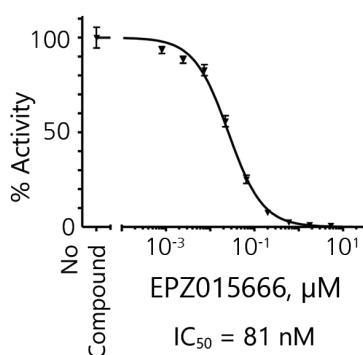
	AlphaLISA™	Chemiluminescent	Colorimetric	Fluorescence Polarization	Fluorogenic	Luminescent	TR-FRET	UV Spectrophotometric	
High specificity	✓	✓	✓	✓	✓	✓	✓	✓	
High sensitivity with low background noise.	✓	✓		✓	✓	✓	✓	✓	
Simple operation	✓		✓	✓	✓	✓	✓		
Homogeneous, no-wash format makes it simple for high throughput screening.	✓		✓	✓	✓	✓	✓		
Real-time readings do not require equilibrium conditions.	✓		✓	✓		✓	✓		
Small volume reactions save on sample consumption.	✓		✓			✓	✓		
Large dynamic range, up to 4-5 logs.	✓	✓			✓				
Highly stable signal allows for repetitive measurements.			✓			✓			
Common, low-cost instrumentation		✓				✓			
Use of lanthanide donor fluorophores minimizes photobleaching.					✓				

Comparison of PRMT5 activity inhibition by EPZ015666 across three assay formats:

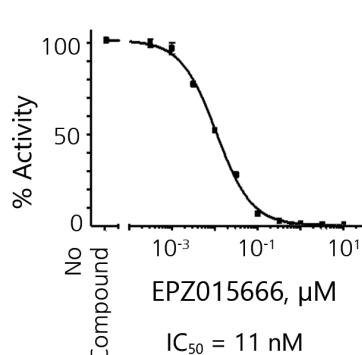
Chemiluminescent Assay



TR-FRET Assay



AlphaLISA® Assay



# Screening & Profiling

BPS Bioscience offers rapid and dependable screening services to determine potency and selectivity of small molecules, peptides, and antibodies against 110+ kinase targets. Our team of experts along with our broad services portfolio make it easy to:

- Screen for inhibitors/targets across full protein families
- Select from IC<sub>50</sub> determination and single point concentrations
- Perform follow-up studies using the same proteins manufactured in-house
- Get questions answered or project guidance in a time-efficient manner

## Our Deliverables to You



### Detailed Results

Extensive report with raw and analyzed data, graphs, and detailed protocols. Typically includes positive controls for inhibition.



### Fast Turnaround

Receive results within weeks after our receipt of your compounds.



### Quality & Consistency

Proteins and enzymes synthesized and QC'd in-house to ensure the highest level of inter- and intra-assay consistency.



### Extensive Selection

Choose from over 200 in-house developed epigenetics-relevant assays, including many novel assay kits.



### Platform Options

Choose the format that works best for your experiment or perform orthogonal testing on multiple platforms to verify your results.

## Key Targets

Deacetylases		Methyltransferases						Demethylases	
HDACs	Sirtuins	DNMTs	EZHs	PRMTs	SETDs	SMYDs	KDMs	TETs	
HDAC1	SIRT1	DNMT1	EZH1	PRMT1	SETDB1	SMYD2	LSD1, LSD2	TET1	
HDAC2	SIRT2	DNMT3A	EZH2	PRMT3	SETD2	SMYD3	FBXL10, FBXL11	TET2	
HDAC3	SIRT3	DNMT3B	EZH2 (A677G)	PRMT4	SET7/9	SMYD4	JMJD1A, JMJD1B		
HDAC4	SIRT5	DNMT3L	EZH2 (A738T)	PRMT5	SET8		JMJD2A, JMJD2B,		
HDAC5	SIRT6		EZH2 del153Y	PRMT6			JMJD2C, JMJD2D,		
HDAC6			EZH2 (P132S)	PRMT7			JMJD2E		
HDAC7			EZH2 (R679C)	PRMT8			JARID1A, JARID1B,		
HDAC8			EZH2 (Y641C)	PRMT9			JARID1C		
HDAC9			EZH2 (Y641F)				UTX, JMJD3		
HDAC10			EZH2 (Y641H)				JHDM1D, PHF8		
HDAC11			EZH2 (Y641N)						
			EZH2 (Y641S)						

Acetyltransferases	Catalog #	Bromodomains	Catalog #
10x HAT Assay Buffer	50095	ATAD2B Inhibitor Screening Assay Kit	32605
AICAR	27049	ATAD2B TR-FRET Assay Kit	32620
ATAT1, GST-tag Recombinant	50072	BAY-299	82063
ATF2, GST-Tag Recombinant	40520	BAZ1A, GST-tag Recombinant	31136
C646	27613-1	BAZ1B, GST-tag Recombinant	31137
C646	27613-2	BAZ1B, His-tag Recombinant	31145
GCN5 (727-837), His-tag Recombinant	31114	BAZ2B Inhibitor Screening Assay Kit	32600
GCN5 (KAT2A), FLAG-tag Recombinant	50074	BAZ2B, His-Tag Recombinant	31113
GCN5 Chemiluminescent Assay Kit	50079L	BET Bromodomain Ligand	33000
HAT Stop Solution	50096	BI-2536	82062
KAT5 (TIP60), GST-Tag Recombinant	79422	BI-9564	82064
p300 (KAT3B), His-GST-tag Recombinant	50071	BPTF/FALZ TR-FRET Assay Kit	32632
P300 Chemiluminescent Assay Kit	79705	BRD1 TR-FRET Assay Kit	32625
P300 Homogeneous Assay Kit	50078	BRD1, GST-tag Recombinant	31011
PCAF (KAT2B), His-tag Recombinant	31120	BRD1, His-tag Recombinant	31010
		BRD2 (BD1), His-tag Recombinant	31022
Bromodomains	Catalog #		
(+)-JQ1	27401	BRD2 (BD1+BD2) TR-FRET Assay Kit	32614
(+)-JQ1	27400	BRD2 (BD1+BD2), GST-tag Recombinant	31024
(+)-JQ1	27402	BRD2 (BD1+BD2), His-tag Recombinant	31025
(-)-JQ1	27335	BRD2 (BD2), GST-tag Recombinant	31023
(-)-JQ1	27334	BRD2 (BD2), His-tag Recombinant	31020
3x BRD Homogeneous Assay Buffer 1	33001	BRD2, GST-tag Recombinant	31021
3x BRD Homogeneous Assay Buffer 2	33007	BRD3 (BD1), GST-tag Recombinant	31032
3x BRD Homogeneous Detection Buffer 2	33006	BRD3 (BD1), His-tag Recombinant	31030
3x BRD TR-FRET Assay Buffer 1	33012	BRD3 (BD1+BD2) TR-FRET Assay Kit	32615
3x BRD TR-FRET Assay Buffer 2	33013	BRD3 (BD1+BD2), GST-tag Recombinant	31035
3x Homogeneous BRD Detection Buffer 1	33002	BRD3 (BD1+BD2), His-tag Recombinant	31034
Anti-BRD2 polyclonal antibody	25211	BRD3 (BD2) Inhibitor Screening Assay Kit	32523
ASH1L, His-tag Recombinant	31147	BRD3 (BD2), GST-tag Recombinant	31033
ATAD2A Inhibitor Screening Assay Kit	32601	BRD3 (BD2), His-tag Recombinant	31031
ATAD2A TR-FRET Assay Kit	32618	BRD4 (BD1) Inhibitor Screening Assay Kit	32514
ATAD2A, GST-tag Recombinant	31121	BRD4 (BD1) TR-FRET Assay Kit	32613
ATAD2A, His-tag Recombinant	31109	BRD4 (BD1), GST-tag Recombinant	31040
ATAD2B (KIAA1240), GST-tag Recombinant	31125	BRD4 (BD1), His-tag Recombinant	31042
ATAD2B (KIAA1240), His-tag Recombinant	31117	BRD4 (BD1+BD2) Recombinant	31047

Bromodomains	Catalog #	Bromodomains	Catalog #
BRD4 (BD1+BD2) TR-FRET Assay Kit	32612	FALZ (BPTF), GST-tag Recombinant	31134
BRD4 (BD1+BD2), GST-Tag Recombinant	31044	FALZ (BPTF), His-tag Recombinant	31131
BRD4 (BD1+BD2), His-Tag Recombinant	31045	GSK-5959	82066
BRD4 (BD2) Inhibitor Screening Assay Kit	32524	GSK-6853	82067
BRD4 (BD2) TR-FRET Assay Kit	32617	I-BET-762	27630-1
BRD4 (BD2), GST-tag Recombinant	31041	I-BET-762	27630-2
BRD4 (BD2), His-tag Recombinant	31043	I-BET151 (GSK1210151A)	27629-1
BRD9 TR-FRET Assay Kit	32621	I-BET151 (GSK1210151A)	27629-2
BRD9, GST-tag Recombinant	31091	I-BET151 (GSK1210151A)	27629-3
BRD9, His-Tag Recombinant	31090	I-BRD9	82068
BRDT (BD1), GST-tag Recombinant	31108	I-CBP112 (Hydrochloride)	27336
BRDT (BD1), His-tag Recombinant	31101	MS-436	82061
BRDT (BD1+BD2) TR-FRET Assay Kit	32616	OTX-015	27636-1
BRDT (BD1+BD2), GST-tag Recombinant	31061	OTX-015	27636-2
BRDT (BD2), GST-tag Recombinant	31060	OTX-015	27636-3
BRDT (BD2), His-tag Recombinant	31100	P300 (1046-1163), His-tag Recombinant	31118
Bromodomain Ligand 2	33003	PB1 (BD4), His-tag Recombinant	31122
Bromodomain Ligand 3	33008	PB1 (BD6), His-tag Recombinant	31133
Bromodomain Ligand 4	33010	PCAF (KAT2B), His-tag Recombinant	31120
Bromodomain Non-Acetylated Ligand 1	33005	PFI-1	27332
Bromodomain Non-Acetylated Ligand 2	33004	PFI-3	82070
Bromodomain Non-Acetylated Ligand 3	33009	PHIP (BD1), GST-tag Recombinant	31142
Bromodomain Non-Acetylated Ligand 4	33011	PHIP (BD1), His-tag Recombinant	31140
Bromosporine	27612	RediSolution™ (+)-JQ1	27403
BRPF1, His-tag Recombinant	31112	RVX-208	27111
BRPF3 TR-FRET Assay Kit	32626	SGC-CBP30	27650-1
BRPF3, GST-tag Recombinant	31130	SGC-CBP30	27650-2
CECR2 Inhibitor Screening Assay Kit	32611	SMARCA2 TR-FRET Assay Kit	40342
CECR2 TR-FRET Assay Kit	32622	SMARCA2, GST-tag Recombinant	31141
CECR2, GST-Tag Recombinant	31138	SMARCA2, His-tag Recombinant	31111
CECR2, His-tag Recombinant	31046	SMARCA2, His-Tag, Biotin-Labeled Recombinant	31129
CHS-828	27333	SMARCA4 (BRG1), GST-tag Recombinant	31132
CREBBP (KAT3A), GST-tag Recombinant	31128	SMARCA4 (BRG1), His-tag Recombinant	31102
CREBBP (KAT3A), His-tag Recombinant	31119	SMARCA4/BRG1 TR-FRET Assay Kit	40341
CREBBP Inhibitor Screening Assay Kit	32607	TAF1 (BD1), His-tag Recombinant	31123
CREBBP TR-FRET Assay Kit	32619	TAF1 (BD1+BD2) TR-FRET Assay Kit	32628

Bromodomains	Catalog #	Deacetylases	Catalog #
TAF1 (BD1+BD2), GST-tag Recombinant	31124	Belinostat (PXD101)	27608-1
TAF1 (BD1/BD2) Inhibitor Screening Assay Kit	32604	Belinostat (PXD101)	27608-2
TAF1 (BD2) Inhibitor Screening Assay Kit	32624	Belinostat (PXD101)	27608-3
TAF1 (BD2) TR-FRET Assay Kit	32627	CBHA	27204
TAF1 (BD2), GST-tag Recombinant	31126	Chidamide	27202
TAF1 (BD2), His-tag Recombinant	31110	CI-994	27021
TAF1L (BD1), GST-tag Recombinant	31105	CUDC-907	27616-1
TAF1L (BD1), His-tag Recombinant	31103	CUDC-907	27616-2
TAF1L (BD1+BD2) TR-FRET Assay Kit	32629	CUDC-907	27616-3
TAF1L (BD1+BD2), GST-tag Recombinant	31107	Droxinostat	27620-1
TAF1L (BD1/BD2) Inhibitor Screening Assay Kit	32603	Droxinostat	27620-2
TAF1L (BD2) Inhibitor Screening Assay Kit	32602	EX527	27103
TAF1L (BD2), GST-tag Recombinant	31106	Fluorogenic HDAC class 2a substrate	50040
TAF1L (BD2), His-tag Recombinant	31104	Fluorogenic HDAC class 2a substrate 2 (green)	50042
TRIM24 (TIF1), GST-tag Recombinant	31127	Fluorogenic HDAC Substrate 2 (Green fluorescence)	50038
TRIM24 (TIF1), His-tag Recombinant	31116	Fluorogenic HDAC substrate 3	50037
TRIM28, GST-tag Recombinant	31139	Fluorogenic HDAC/SIRT substrate 1	50032
TRIM28, His-Tag Recombinant	31146	Fluorogenic Rat HDAC6 Assay Kit	78849
WDR9/BRWD1 (BD2), GST-tag Recombinant	31135	Fluorogenic Sirtuin 5 Substrate	50126
WDR9/BRWD1 (BD2), His-tag Recombinant	31115	Fluorogenic Sirtuin 6 Substrate	50866
WDR9/BRWD1 TR-FRET Assay Kit	32631	HC Toxin	27209
ZMYND11, GST-tag Recombinant	31144	HDAC Assay Buffer	50031
ZMYND8, GST-tag Recombinant	31143	HDAC developer (2X)	50030
		HDAC Fluorogenic Assay Kit (Green)	50034
Deacetylases	Catalog #		
4-iodo-SAHA	27205	HDAC1 Fluorogenic Assay Kit	50061
AGK2	27228	HDAC1 Kinetic Assay Kit	53001
Anti-HDAC1 monoclonal antibody	25286	HDAC1, FLAG-Tag, His-Tag Recombinant	50051
Anti-HDAC1 polyclonal antibody	25287	HDAC1, His-Tag, FLAG-Tag (Mouse) Recombinant	50058
Anti-HDAC2 monoclonal antibody	25288	HDAC10, FLAG-Tag Recombinant	50060
Anti-HDAC3 monoclonal antibody	25289	HDAC11 Fluorogenic Assay Kit	50687
Anti-SIRT1 monoclonal antibody	25316	HDAC11 Recombinant	50021
Apicidin	27329	HDAC2 Fluorogenic Assay Kit	50062
AR-42 (OSU-HDAC42)	27603-1	HDAC2 Kinetic Assay Kit	53002
AR-42 (OSU-HDAC42)	27603-2	HDAC2, FLAG-tag Recombinant	50052
AR-42 (OSU-HDAC42)	27603-3	HDAC2, His-tag Recombinant	50002
		HDAC3 Fluorogenic Assay Kit	50073

Deacetylases	Catalog #	Deacetylases	Catalog #
HDAC3 Kinetic Assay Kit	53003	Mocetinostat	27019
HDAC3/NcoR2, His-Tag, GST-Tag Recombinant	50003	Mocetinostat (MGCD0103)	27109
HDAC4 Fluorogenic Assay Kit	50064	MS-275 (Entinostat)	27011
HDAC4, GST-Tag Recombinant	100339	NcoR2 (SMRT), GST-tag Recombinant	50020
HDAC4, GST-tag, His-tag Recombinant	50004	Oxamflatin	27326
HDAC5 (full length), GST-tag Recombinant	50045	Oxamflatin	27201
HDAC5 Fluorogenic Assay Kit	50065	Panobinostat	27020
HDAC5, Catalytic Domain, His-tag Recombinant	50005	Parthenolide	27768-1
HDAC5, GST-Tag, His-Tag Recombinant	100102	Parthenolide	27768-2
HDAC6 (H216A), GST-Tag Recombinant	50046	PCI 34051	27316
HDAC6 (H611A), GST-tag Recombinant	50066	PCI 34051	27317
HDAC6 Fluorogenic Assay Kit	50076	PCI-24781 (CRA-024781)	27769-1
HDAC6, FLAG-Tag Recombinant	50056	PCI-24781 (CRA-024781)	27769-2
HDAC6, GST-Tag (HEK293-derived) Recombinant	100051	PCI-24781 (CRA-024781)	27769-3
HDAC6, GST-Tag (Mouse) Recombinant	50057	pf-HDAC1, Malaria, His-tag, FLAG-tag Recombinant	50063
HDAC6, GST-Tag (Rat) Recombinant	101332	Pracinostat (SB939)	27771-1
HDAC6, GST-tag (Sf9-derived) Recombinant	50006	Pracinostat (SB939)	27771-2
HDAC7 Fluorogenic Assay Kit	50067	Pracinostat (SB939)	27771-3
HDAC7, GST-tag Recombinant	50007	Pyroxamide	27328
HDAC8 Fluorogenic Assay Kit	50068	Resveratrol	27642-3
HDAC8, His-Tag Recombinant	50008	RGFP966	27643-1
HDAC9 Fluorogenic Assay Kit	50069	RGFP966	27643-2
HDAC9, His-tag Recombinant	50009	RGFP966	27643-3
HNHA	27206	Rocilinostat (ACY-1215)	27644-1
ITF 2357	27327	Rocilinostat (ACY-1215)	27644-2
JNJ-26481585	27632-1	Rocilinostat (ACY-1215)	27644-3
JNJ-26481585	27632-2	Romidepsin (FK228, depsipeptide)	27646-1
JNJ-26481585	27632-3	Romidepsin (FK228, depsipeptide)	27646-2
Kaempferol	27757	SAHA (Vorinostat)	27006
Kinetic HDAC Assay Developer	53029	Salermide	27324
LAQ824 (NVP-LAQ824,Dacinostat)	27760-1	SB 939	27331
LAQ824 (NVP-LAQ824,Dacinostat)	27760-2	Scriptaid	27200
LpxC (E. coli) Recombinant	11128	SIRT1 (Sirtuin1) Fluorogenic Assay Kit	50081
M344	27203	SIRT1/2 Inhibitor IV	27323
MC1568	27761-1	SIRT2 (Sirtuin2) Fluorogenic Assay Kit	50087
MC1568	27761-2	SIRT3 (Sirtuin3) Fluorogenic Assay Kit	50088

Deacetylases	Catalog #	Demethylases	Catalog #
SIRT5 (Sirtuin5) Fluorogenic Assay Kit	50085	4x HDM Assay Buffer 1	52406
Sirtinol	27652-1	4x HDM Assay Buffer 2	52407
Sirtinol	27652-2	4x HDM Assay Buffer 3	52408
Sirtuin 1 (193-741), GST-tag Recombinant	50012	4x HDM Assay Buffer 4	52409
Sirtuin 2, His-Tag Recombinant	50013	4x HDM Incomplete Assay Buffer	52410
Sirtuin 3, GST-Tag Recombinant	50014	ALKBH5 Chemiluminescent Assay Kit	79659
Sirtuin 4, GST-tag Recombinant	50015	ALKBH5, FLAG-Tag Recombinant	100057
Sirtuin 5, GST-tag Recombinant	50016	Anti-Jarid1b polyclonal antibody	25293
Sirtuin 6 Fluorogenic Assay Kit	50022	Anti-Jarid1c polyclonal antibody	25294
Sirtuin 6, GST-tag Recombinant	50017	Anti-JARID2 polyclonal antibody	25295
Sirtuin 7, FLAG-Tag Recombinant	50018	Anti-JMJD2a polyclonal antibody	25296
Sirtuin Assay Buffer	50090	Anti-JMJD2b polyclonal antibody	25297
Sirtuin Developer (2X)	50089	Anti-JMJD2D (KDM4D) polyclonal antibody	25300
Sodium Phenylbutyrate	27654	Anti-LSD1 polyclonal antibody	25302
Splitomicin	27330	Anti-LSD2 (KDM1B) polyclonal antibody	25299
SRT1720	27655-1	Anti-PHF8 polyclonal antibody	25308
SRT1720	27655-2	Anti-TET2 monoclonal antibody	25320
SRT1720	27655-3	Biotinylated Histone H3 Peptide Substrate	79656
Suberohydroxamic Acid (SBHA)	27210	Blocking Buffer 1 (Multiple Epigenetics Assay Kits)	79556
Tenovin-1	27660-1	Blocking Buffer 4 (Multiple Epigenetics Assay Kits)	52100
Tenovin-1	27660-2	FBXL10 (KDM2B) Chemiluminescent Assay Kit	50157
Tenovin-1	27660-3	FBXL10 (KDM2B) Homogeneous Assay Kit	50610
Tenovin-6	27661-1	FBXL10 (KDM2B, JHDM1B), FLAG-tag Recombinant	50120
Tenovin-6	27661-2	FBXL11 (KDM2A) Homogeneous Assay Kit	50611
Tenovin-6	27661-3	FBXL11 (KDM2A, JHDM1A), FLAG-Avi-tags Recombinant	50156
Trichostatin A (TSA)	27102	FTO Chemiluminescent Assay Kit	79344
Tubacin	27662	FTO, His-Tag (Sf9 derived) Recombinant	79480
Tubastatin A	27663-1	FTO, His-Tag Recombinant	79306
Tubastatin A	27663-2	GSK J4 HCl	27627-1
Tubastatin A	27663-3	GSK J4 HCl	27627-2
Tubastatin A hydrochloride	27108	H3K4me1 Substrate	50101
Valproic Acid	27207	H3K4me2 Substrate	50158
<b>Demethylases</b>		<b>Catalog #</b>	
JMJD2D (KDM4D) Chemiluminescent Assay Kit	50418	Histone H3 K27cMe1, Full Length, His-tag Recombinant	52081
4x Detection Buffer 1	52301	Histone H3 K27cMe2, Full Length, His-tag Recombinant	52082
		Histone H3 K27cMe3, Full Length, His-tag Recombinant	52083
		Histone H3 peptide (1-21), Dimethyl K9, Biotin-labeled Recombinant	50351

Demethylases	Catalog #	Demethylases	Catalog #
Histone H3 peptide (1-21), Monomethyl K9, Biotin-labeled Recombinant	50353	JMJD3 (KDM6B) Chemiluminescent Assay Kit	50406
Histone H3 peptide (1-21), Trimethyl K9, Biotin-labeled Recombinant	50350	JMJD3 (KDM6B) Homogeneous Assay Kit	50416
Histone H3-H4 Tetramer H3K27cMe2, His-tag Recombinant	52086	JMJD3 (KDM6B), FLAG-tag Recombinant	50115
Histone H3-H4 Tetramer H3k27cMe3, His-tag Recombinant	52087	JMJD6, GST-tag Recombinant	50140
Histone Octamer H3K27cMe2, His-tag Recombinant	52090	LSD1 (KDM1A) Chemiluminescent Assay Kit	50109
Histone Octamer H3K27cMe3, His-tag Recombinant	52091	LSD1 (KDM1A) Homogeneous Assay Kit	50108
JARID1A (KDM5A) Chemiluminescent Assay Kit	50513	LSD1 (KDM1A), Variant 2 Recombinant	50097
JARID1A (KDM5A) Homogeneous Assay Kit	50510	LSD1 (KDM1A), Variant 2, GST-tag Recombinant	50100
JARID1A (KDM5A), FLAG-tag Recombinant	50110	LSD2 (KDM1B), GST-Tev-Tag Recombinant	50124
JARID1A (KDM5A), His-Avi-tags, FLAG-tag Recombinant	50155	LSD2(KDM1B) Homogeneous Assay Kit (Dimethyl)	50614
JARID1B (KDM5B) Chemiluminescent Assay Kit	50517	LSD2(KDM1B) Homogeneous Assay Kit (Monomethyl)	50613
JARID1B (KDM5B) Homogeneous Assay Kit	50512	Mouse JARID1B (KDM5B, PLU-1), FLAG-tag Recombinant	50122
JARID1B (KDM5B) TR-FRET Assay Kit	50514	OG-L002	27635-1
JARID1B (KDM5B), Full Length, FLAG-tag Recombinant	50128	OG-L002	27635-2
JARID1B (KDM5B, PLU-1), His-FLAG-tags Recombinant	50121	PHF8 (KDM7B) Homogeneous Assay Kit	50515
JARID1C (KDM5C) Homogeneous Assay Kit	50511	PHF8 (KDM7B), (2-525), FLAG-tag Recombinant	50131
JARID1C(KDM5C), FLAG-tag Recombinant	50112	PHF8 (KDM7B), Full-length, FLAG-tag Recombinant	50132
JHDM1D (KDM7A) Chemiluminescent Assay Kit	50612	Posaconazole	27639-1
JHDM1D (KDM7A) Homogeneous Assay Kit	50420	Posaconazole	27639-2
JHDM1D (KDM7A), FLAG-tag Recombinant	50419	Secondary Anti-HRP-labeled Antibody 2	52131H
JMJD1A (KDM3A), FLAG-tag Recombinant	50130	TET1 Chemiluminescent Assay Kit	50651
JMJD1A Homogeneous Assay Kit	50412	TET1, FLAG-tag Recombinant	50161
JMJD1B (KDM3B), FLAG-tag Recombinant	50421	TET2 Chemiluminescent Assay Kit	50652
JMJD1B Homogeneous Assay Kit	79839	TET2, FLAG-Tag, His-Tag (916-2002) Recombinant	78559
JMJD2A (KDM4A) (1-350), His-Tag Recombinant	50123	TET2, FLAG-Tag, His-Tag (catalytic domain) Recombinant	100408
JMJD2A (KDM4A) Homogeneous Assay Kit	50413	TET3, FLAG-Tag, His-Tag Recombinant	50163
JMJD2B (KDM4B) Homogenous Assay Kit	50414	Tranylcypromine (2-PCPA)	27305
JMJD2B (KDM4B), GST-tag Recombinant	50111	UTX (KDM6A), C-term FLAG-tag Recombinant	50116
JMJD2B(KDM4B) , FLAG-tag Recombinant	50104	UTX (KDM6A), N-terminal FLAG-tag Recombinant	50119
JMJD2C (KDM4C) Chemiluminescent Assay Kit	50405	UTX Chemiluminescent Assay Kit	50615
JMJD2C (KDM4C) Homogeneous Assay Kit	50415	UTX(KDM6A) Homogeneous Assay Kit	50516
JMJD2C, GST-Tag Recombinant	50105		
JMJD2D (KDM4D), GST-tag Recombinant	50117	Histones	Catalog #
JMJD2D Homogeneous Assay Kit	79838	Anti-H1.0 polyclonal antibody	25225
JMJD2E (KDM4DL) Homogeneous Assay Kit	50417	Anti-H2A pan polyclonal antibody	25226
JMJD2E (KDM4DL), GST-tag Recombinant	50118		

Histones	Catalog #	Histones	Catalog #
Anti-H2A.XS139p polyclonal antibody	25227	Anti-H3K9ac monoclonal antibody	25265
Anti-H2A.Zac polyclonal antibody	25228	Anti-H3K9acS10p polyclonal antibody	25267
Anti-H2AK119ub polyclonal antibody	25229	Anti-H3K9me1 polyclonal antibody	25268
Anti-H2AK5ac polyclonal antibody	25230	Anti-H3K9me2 monoclonal antibody	25269
Anti-H2B pan polyclonal antibody	25231	Anti-H3K9me2 polyclonal antibody	25270
Anti-H2BK12ac polyclonal antibody	25232	Anti-H3K9me3 monoclonal antibody	25271
Anti-H2BK15ac polyclonal antibody	25233	Anti-H3K9me3 polyclonal antibody	25272
Anti-H2Bs14p monoclonal antibody	25234	Anti-H3K9me3S10p polyclonal antibody	25273
Anti-H3 pan polyclonal antibody	25235	Anti-H3K9un monoclonal antibody	25274
Anti-H3K18ac polyclonal antibody	25236	Anti-H3R17me2 polyclonal antibody	25275
Anti-H3K27ac monoclonal antibody	25238	Anti-H3R17me2(asym)K18ac polyclonal antibody	25276
Anti-H3K27ac polyclonal antibody	25239	Anti-H3S10p polyclonal antibody	25277
Anti-H3K27me1 polyclonal antibody	25240	Anti-H4 pan polyclonal antibody	25278
Anti-H3K27me2 polyclonal antibody	25241	Anti-H4K20me1 monoclonal antibody	25279
Anti-H3K27me3 monoclonal antibody	25243	Anti-H4K20me1 polyclonal antibody	25280
Anti-H3K27me3 polyclonal antibody	25244	Anti-H4K20me3 polyclonal antibody	25281
Anti-H3K27me3S28p polyclonal antibody	25245	Anti-H4K5,8,12,16ac polyclonal antibody	25282
Anti-H3K36me1 polyclonal antibody	25246	Anti-H4K5,8,12ac polyclonal antibody	25283
Anti-H3K36me2 monoclonal antibody	25247	Anti-H4K5ac polyclonal antibody	25284
Anti-H3K36me2 polyclonal antibody	25248	Anti-H4K8ac polyclonal antibody	25285
Anti-H3K36me3 monoclonal antibody	25249	Anti-H4R3me2(sym) polyclonal antibody	78558
Anti-H3K36me3 polyclonal antibody	25250	Histone H3 peptide (1-21), Dimethyl K9, Biotin-labeled Recombinant	50351
Anti-H3K4ac polyclonal antibody	25251	Histone H4 Peptide (1-23), Monomethyl R3, Biotin-labeled Recombinant	51201
Anti-H3K4me1 monoclonal antibody	25252	Histone Octamer H3K27cMe2, His-tag Recombinant	52090
Anti-H3K4me1 polyclonal antibody	25253		
Anti-H3K4me2 monoclonal antibody	25254		
Anti-H3K4me2 polyclonal antibody	25255	<b>Methyl-lysine Reader(s)</b>	<b>Catalog #</b>
Anti-H3K4me3 monoclonal antibody	25256	Anti-CBX2 polyclonal antibody	25212
Anti-H3K4me3 polyclonal antibody	25257	Anti-Chd1 polyclonal antibody	25213
Anti-H3K4un monoclonal antibody	25258	CBX1 Inhibitor Screening Assay Kit	79052
Anti-H3K56ac polyclonal antibody	25259	CBX1, GST-tag Recombinant	55009
Anti-H3K64me3 polyclonal antibody	25260	CBX2, GST-tag Recombinant	55011
Anti-H3K79me1 polyclonal antibody	25261	CBX2, His-tag Recombinant	55012
Anti-H3K79me2 polyclonal antibody	25262	CBX7, His-tag Recombinant	55017
Anti-H3K79me3 polyclonal antibody	25263	CDY1, GST-tag Recombinant	55007
Anti-H3K9/14ac polyclonal antibody	25264	CDY1, His-tag Recombinant	55015
		CHD2, GST-tag Recombinant	55005

Methyl-lysine Reader(s)	Catalog #	Methyltransferases	Catalog #
L3MBTL1 Inhibitor Screening Assay Kit	55100	Anti-DNMT1 polyclonal antibody	25216
L3MBTL1 TR-FRET Assay Kit	55200	Anti-DNMT2 polyclonal antibody	25217
L3MBTL1, GST-tag Recombinant	55000	Anti-DNMT3A polyclonal antibody	25218
L3MBTL1, His-tag Recombinant	55002	Anti-DNMT3B polyclonal antibody	25219
L3MBTL2, His-tag Recombinant	55018	Anti-EHMT1 polyclonal antibody	25220
L3MBTL3, GST-tag Recombinant	55013	Anti-EZH2 monoclonal antibody	25222
L3MBTL3, His-tag Recombinant	55014	Anti-EZH2 polyclonal antibody	25223
UHRF1 (108-286), GST-tag Recombinant	55003	Anti-G9a polyclonal antibody	25224
UHRF1 (108-286), His-tag Recombinant	55004	Anti-RbAp 46/48 monoclonal antibody	25311
UHRF1, His-FLAG-Tags Recombinant	55001	Anti-Set9 polyclonal antibody	25312
UNC1215	27406	Anti-Setd1a polyclonal antibody	25313
UNC1215	27405	Anti-Setd1b polyclonal antibody	25314
UNC669	27664-1	Anti-SETD8 polyclonal antibody	25315
UNC669	27664-2	Anti-WDR5 polyclonal antibody Anti- Adenosyl Homocysteine Antibody	25321 52030
Methyltransferases	Catalog #		
3-Deazaneplanocin (DZNep)	27617-1	Ash2L, His-tag Recombinant	71201
3-Deazaneplanocin (DZNep)	27617-2	Azacytidine (Vidaza)	27107
3-Deazaneplanocin A (DZNep) hydrochloride	27618-1	BIX01294 (hydrochloride hydrate)	27220
3-Deazaneplanocin A (DZNep) hydrochloride	27618-2	Blocking Buffer 1 (Multiple Epigenetics Assay Kits)	79556
400 µM S-Adenosylmethionine	52120	Blocking Buffer 4 (Multiple Epigenetics Assay Kits)	52100
4x Detection Buffer	79389	Chaetocin	27221
4x DNMT Assay Buffer 1	52200	CoREST, His-tag Recombinant	50274
4x DNMT Assay Buffer 2	52201	CTPB	27050
4x HMT Assay Buffer 1	52160	DNMT Universal Chemiluminescent Assay Kit	52035
4x HMT Assay Buffer 1	52160-A	DNMT1 Chemiluminescent Assay Kit	52050L
4x HMT Assay Buffer 2	52170	DNMT1, GST-tag Recombinant	51101
4x HMT Assay Buffer 2A	52170-A	DNMT1, His-FLAG-tag Recombinant	51100
4x HMT Assay Buffer 3	52180	DNMT2, GST-tag Recombinant	51102
4x HMT Assay Buffer 4	52190	DNMT3A Chemiluminescent Assay Kit	52033
4x HMT Assay Buffer 5	52191	DNMT3A, GST-Tag Recombinant	51103
4x HMT Assay Buffer 6	52192	DNMT3B Chemiluminescent Assay Kit	52034
4x HMT Assay Buffer 7	52193	DNMT3B, His-tag, FLAG-tag (Mouse) Recombinant	51105
4x METTL Assay Buffer	78368	DNMT3B/DNMT3L Complex, GST-tag, His-tag Recombinant	51109
AEBP2, His-tag Recombinant	50281	DNMT3B/DNMT3L Complex, His-tag, GST-tag Recombinant	51104
AMI-1 (sodium salt)	27307	DOT1L (KMT4) Chemiluminescent Assay Kit	52202

Methyltransferases	Catalog #	Methyltransferases	Catalog #
DOT1L (KMT4), GST-tag Recombinant	51005	EZH2 (R679C)/EED/SUZH12/RbAp48/AEBP2, FLAG-tag, His-tag Recombinant	51033
DPY30, His-tag Recombinant	71203	EZH2 (Y641C) /EED/SUZ12/RbAp48/AEBP2, His-tag, FLAG-tag Recombinant	51029
EED, FLAG-tag Recombinant	50280	EZH2 (Y641F) /EED/SUZ12/RbAp48/AEBP2, His-tag, FLAG-tag Recombinant	51017
EHMT1 (GLP) Chemiluminescent Assay Kit	53007	EZH2 (Y641F) Chemiluminescent Assay Kit	52075
EHMT1 (GLP), GST-tag Recombinant	51020	EZH2 (Y641F) Chemiluminescent Assay Kit	79276
Ellagic acid	27219	EZH2 (Y641H) /EED/SUZ12/RbAp48/AEBP2, His-tag, FLAG-tag Recombinant	51011
Entacapone	27621-1	EZH2 (Y641N) /EED/SUZ12/RbAp48/AEBP2, His-Tag, FLAG-Tag Recombinant	51028
Entacapone	27621-2	EZH2 (Y641N) Chemiluminescent Assay Kit	79277
EPZ-6438	27626-1	EZH2 (Y641N) TR-FRET Assay Kit	52076
EPZ-6438	27626-2	EZH2 (Y641S) /EED/SUZ12/RbAp48/AEBP2, His-tag, FLAG-tag Recombinant	51013
EPZ-6438	27626-3	EZH2 / EED / SUZ12, His-Tag, FLAG-Tag Recombinant	51003
EPZ004777	27622	EZH2 Chemiluminescent Assay Kit	52009L
EPZ005687	27624-1	EZH2(Y641F) TR-FRET Assay Kit	52085
EPZ005687	27624-2	EZH2-EED Binding Assay Kit	52059
EPZ015666	79043-1	EZH2/EED, His-tag, FLAG-tag Recombinant	52066
EPZ015666	79043-2	EZH2/EED/SUZ12/RbAp48, His-tag, FLAG-tag Recombinant	51002
EPZ5676	27625-1	EZH2/EED/SUZ12/RbAp48/AEBP2, FLAG-tag, His-tag Recombinant	51099
EPZ5676	27625-2	EZH2/Homogeneous Assay Kit	51004
EZH1 Chemiluminescent Assay Kit	52079	EZH2/EED/SUZ12/RbAp48/AEBP2, FLAG-tag, His-tag Recombinant	51098
EZH1 Chemiluminescent Assay Kit	52990	EZH2/SUZ12/RbAp48, His-GST-tags, His-tag Recombinant	51000
EZH1/EED/SUZ12, His-tag, FLAG-tag Recombinant	51006	G9a (EHMT2), GST-tag (E. coli-derived) Recombinant	51001
EZH1/EED/SUZ12/RbAp48/AEBP2, His-tag, FLAG-tag Recombinant	51007	G9a (EHMT2), GST-tag (Sf9-derived) Recombinant	52001L
EZH2 (A677G) Chemiluminescent Assay Kit	52077	G9a Chemiluminescent Assay Kit	52051
EZH2 (A677G) Chemiluminescent Assay Kit	52070	G9a Homogeneous Assay Kit	52043
EZH2 (A677G)/EED/SUZ12/RbAp48/AEBP2, His-tag, FLAG-tag Recombinant	51037	Histone H1, Full Length, His-tag Recombinant	52049
EZH2 (A677T)/EED/SUZ12/RbAp48/AEBP2, His-tag, FLAG-tag Recombinant	51039	Histone H2a, Full Length, Biotin-labeled, His-tag Recombinant	52036
EZH2 (A687V)/EED/SUZ12/RbAp48/AEBP2, His-tag, FLAG-tag Recombinant	51038	Histone H2a, Full Length, His-tag Recombinant	52025
EZH2 (A738T) /EED/SUZ12/RbAp48/AEBP2, FLAG-Tag, His-Tag Recombinant	51061	Histone H2a, Full Length, His-tag Recombinant	52021
EZH2 (A738T) Chemiluminescent Assay Kit	52067	Histone H2b, Full Length, His-tag Recombinant	52026
EZH2 (A738T) Chemiluminescent Assay Kit	79275	Histone H2b, Full Length, His-tag Recombinant	52022
EZH2 (del153Y)/EED/SUZ12/RbAp48/AEBP2, FLAG, His-tag Recombinant	51027	Histone H3 (2-58), GST-tag Recombinant	52013
EZH2 (E740K)/EED/SUZH12/RbAp48/AEBP2, FLAG-tag, His-tag Recombinant	51032	Histone H3 (K27M), His-tag Recombinant	52101
EZH2 (F667I)/EED/SUZH12/RbAp48/AEBP2, FLAG-tag, His-tag Recombinant	51035	Histone H3 K27C, Full Length, His-tag Recombinant	52080
EZH2 (H689Y)/EED/SUZ12/RbAp48/AEBP2, FLAG-tag, His-tag Recombinant	51023	Histone H3 K27cMe1, Full Length, His-tag Recombinant	52081
EZH2 (non-complexed), His-GST-tags Recombinant	50279		
EZH2 (P132S)/EED/SUZ12/RbAp48/AEBP2, FLAG-tag, His-tag Recombinant	51019		

Methyltransferases	Catalog #	Methyltransferases	Catalog #
Histone H3 K27cMe2, Full Length, His-tag Recombinant	52082	MLL1 Complex Chemiluminescent Assay Kit	53008
Histone H3 peptide (1-21)	52010	MLL1/WDR5/Ash2L/RbBP5 (MLL1/WAR Complex), His-tag Recombinant	51022
Histone H3 peptide (1-21), Biotin-labeled Recombinant	52011	MLL1/WDR5/Ash2L/RbBP5/DPY30 (MLL1/WARD Complex), FLAG, His-Tag Recombinant	51021
Histone H3 peptide (21-44)	52016	MLL3 Complex Chemiluminescent Assay Kit	79758
Histone H3 peptide (21-44), Biotin-labeled Recombinant	52017	MLL3/WDR5/Ash2L/RbBP5/DPY30 Complex Recombinant	100323
Histone H3 Peptide (21-44), Monomethyl K27, Biotin-labeled Recombinant	52020	MLL4 Complex Chemiluminescent Assay Kit	79757
Histone H3 Recombinant	79055-2	Native Core Histone Octamers	52065
Histone H3 Recombinant	79055-1	Native Nucleosome (HEK293-derived)	52015
Histone H3(K9) Universal Methyltransferase Assay Kit	52072	Native Nucleosome (HeLa-derived)	52039
Histone H3, Full Length, Biotin-labeled, His-tag	52053	NNMT, His-Tag Recombinant	11357
Histone H3, Full Length, Biotin-labeled, His-tag	52047	NSD1 Chemiluminescent Assay Kit	53010
Histone H3, Full Length, His-tag Recombinant	52027	NSD1, GST-tag Recombinant	51024
Histone H3, Full Length, His-tag Recombinant	52023	NSD2 (catalytic domain), GST-tag Recombinant	51026
Histone H3-H4 Tetramer H3K27C Unmodified, His-tag Recombinant	52084	NSD2 Chemiluminescent Assay Kit	53009
Histone H3-H4 Tetramer H3K27cMe2, His-tag Recombinant	52086	NSD2 Chemiluminescent Assay Kit	79359
Histone H3/H4 Tetramer (full length), His-tag Recombinant	52064	NSD3 (1021-1322) (WHSC1L1), GST-tag Recombinant	51036
Histone H4 (2-58), GST-tag Recombinant	52019	NSD3 Chemiluminescent Assay Kit	53012
Histone H4 peptide (1-21)	52014	NSD3 Chemiluminescent Assay Kit	79358
Histone H4 peptide (1-21), Biotin-labeled Recombinant	52018	PCGF6, GST-tag Recombinant	55016
Histone H4 peptide (1-21), Dimethyl R3, Biotin-labeled Recombinant	51200	PRDM12, GST-tag Recombinant	50320
Histone H4(R3) Methyltransferase Assay Kit	52074	PRDM14, GST-tag Recombinant	50321
Histone H4, Full Length, His-tag Recombinant	52024	PRDM9 Chemiluminescent Assay Kit	79625
Histone H4, Full Length, His-tag Recombinant	52028	PRDM9, GST-Tag Recombinant	100078
Histone Octamer (Full length) Recombinant	52040	PRMT1 Chemiluminescent Assay Kit	52004L
Histone Octamer (full length), His-tag Recombinant	52037	PRMT1 Homogeneous Assay Kit	52054
Histone Octamer H3K27C Unmodified, His-tag Recombinant	52088	PRMT1, GST-Tag (E. coli-derived) Recombinant	51040
Histone Octamer H3K27cMe2, His-tag Recombinant	52090	PRMT1, His-GST-tag (Sf9-derived) Recombinant	51041
Histone Octamer, Biotin-labeled, His-tag Recombinant	52061	PRMT2, FLAG-Tag Recombinant	100471
Histone-H2A-(107-122)-Ac-OH	27628-1	PRMT3 Chemiluminescent Assay Kit	52005L
Histone-H2A-(107-122)-Ac-OH	27628-2	PRMT3 Homogeneous Assay Kit	52055
MECP2, GST-tag Recombinant	50250	PRMT3, GST-tag Recombinant	51043
Methnase (SETMAR), FLAG-tag Recombinant	51018	PRMT4 (CARM1), FLAG-tag (HEK293-derived) Recombinant	51047
METTL3, FLAG-Tag Recombinant	100055	PRMT4 Chemiluminescent Assay Kit	52041L
METTL3/METTL14 Complex Chemiluminescent Assay Kit	79614	PRMT4 Homogeneous Assay Kit	52068
METTL3/METTL14 Protein Complex, FLAG-Tag, His-Tag Recombinant	100105	PRMT4 peptide substrate, Biotin-labeled	50352
MGMT, His-Tag Recombinant	101602	PRMT5 Chemiluminescent Assay Kit	52002L

Methyltransferases	Catalog #	Methyltransferases	Catalog #
PRMT5 Chemiluminescent Assay Kit	52073	SMYD2 (KMT3C) Chemiluminescent Assay Kit	51031
PRMT5 Homogeneous Assay Kit	52052	SMYD2 (KMT3C), FLAG-Tag Recombinant	51014
PRMT5 TR-FRET Assay Kit	52171	SMYD3 (35-end), GST-tag Recombinant	51015
PRMT5, FLAG-Avi-Tag, Biotin-Labeled Recombinant	79191	SMYD3 (full length), FLAG-tag Recombinant	51016
PRMT5/MEP50, Biotin-Labeled Recombinant	52102	SMYD3 Homogeneous Assay Kit	52063
PRMT5/MEP50, FLAG-Tag, His-Tag (HEK293-derived) Recombinant	51045	SMYD4 Chemiluminescent Assay Kit	53013
PRMT5/MEP50, FLAG-Tag, His-Tag (Sf9-derived) Recombinant	51048	SMYD4, FLAG-tag Recombinant	51057
PRMT6 (HRMT1L6) Chemiluminescent Assay Kit	52046	SUV39H1 Chemiluminescent Assay Kit	52006L
PRMT6 (HRMT1L6) Homogeneous Assay Kit	52056	SUV39H1, GST-tag Recombinant	51070
PRMT6, His-tag Recombinant	51049	SUV39H2 Activity Assay Kit	52008L
PRMT7 Homogeneous Assay Kit (Automethylation)	79721	SUV39H2, GST-Tag Recombinant	51080
PRMT7 Homogeneous Assay Kit (Histone H2b)	79720	SUV4-20H1 Chemiluminescent Assay Kit	53011
PRMT7, His-FLAG-tag Recombinant	51054	SUV4-20H1, GST-tag Recombinant	51090
PRMT8 Chemiluminescent Assay Kit	52062	SUV4-20H2, GST-tag Recombinant	51060
PRMT8 Homogeneous Assay Kit	52058	SUZ12, His-FLAG-tags Recombinant	50282
PRMT8, His-FLAG-tags Recombinant	51052	UNC0224	27303
PRMT9 Homogeneous Assay Kit	52069	UNC0321 (TFA salt)	27306
PRMT9, FLAG-Tag, His-Tag Recombinant	79124	UNC0638	27222
PRMT9, FLAG-Tag, Strep-Tag Recombinant	52057	WDR5, FLAG-tag Recombinant	51111
RbAp48, His-tag Recombinant	50297	WDR5, His-tag Recombinant	71200
RbBP5, His-tag Recombinant	71202	WDR5, His-Tag, Biotin-Labeled Recombinant	80016
Recombinant Nucleosome (E. coli-derived)	52038	<b>RNA Epigenetics</b>	
Recombinant Nucleosome (E. coli-derived)	52333	ADAR1, FLAG-Tag Recombinant	100472
Recombinant Nucleosome (E. coli-derived), Biotin-labeled, His-Tag	52048	ADAR2 (ADARB1), FLAG-Tag Recombinant	101164
S-adenosyl homocysteine	52031	ALKBH5 Chemiluminescent Assay Kit	79659
S-adenosyl homocysteine, Fluorescein-labeled	52012	ALKBH5, FLAG-Tag Recombinant	100057
SAP145, GST-Tag Recombinant	79111	Anti-DNMT2 polyclonal antibody	25217
Secondary Anti-HRP-labeled Antibody 2	52131H	DNMT2, GST-tag Recombinant	51102
Set7/9 (Sf9), His-GST-tags Recombinant	50019	FTO Chemiluminescent Assay Kit	79344
Set7/Set9, GST-tag (E. coli-derived) Recombinant	51010	FTO, His-Tag (Sf9 derived) Recombinant	79480
Set8 Chemiluminescent Assay Kit	82244	FTO, His-Tag Recombinant	79306
Set8, Active, GST-Tag Recombinant	51008	Lin28A, FLAG-Avi-Tag Recombinant	79254
SETD2 Chemiluminescent Assay Kit	52060	Lin28A, FLAG-Avi-Tag, Biotin-Labeled Recombinant	79072
SETD2, GST-tag Recombinant	53019	Lin28B, FLAG-Avi-Tag Recombinant	79255
SetDB1 Chemiluminescent Assay Kit	51056	Lin28B, FLAG-Avi-Tag, Biotin-Labeled Recombinant	79073
SetDB1, His-GST-tags Recombinant	51055		

RNA Epigenetics	Catalog #
METTL3, FLAG-Tag Recombinant	100055
METTL3/METTL14 Complex Chemiluminescent Assay Kit	79614
METTL3/METTL14 Protein Complex, FLAG-Tag, His-Tag Recombinant	100105
TUT4, FLAG-tag (Human) Recombinant	71259
TUT4, FLAG-tag, Biotin-Labeled (Human) Recombinant	71258



## FIND A TRUSTED PARTNER FOR YOUR NEXT PROJECT



Ordering Information  
[orders@bpsbioscience.com](mailto:orders@bpsbioscience.com)



Technical Support  
[support@bpsbioscience.com](mailto:support@bpsbioscience.com)



US Sales Support  
[sales-team@bpsbioscience.com](mailto:sales-team@bpsbioscience.com)



International Sales Support  
[international@bpsbioscience.com](mailto:international@bpsbioscience.com)

BPS Bioscience, Inc.  
6405 Mira Mesa Blvd, Suite 100  
San Diego, CA 92121  
Tel: 858-202-1401

[bpsbioscience.com](http://bpsbioscience.com)

