

# diagenode

A Hologic Company

## SIMPLY SUPERIOR EPIGENETIC ANTIBODIES







Click on the antibody's name to visit the product page

## Unmodified histones and pan antibodies



PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS
<a href="#">H2A pan pAb</a>	C15410166	50 µg	Hu	ChIP - ELISA - WB - IF
<a href="#">H2B pan pAb</a>	C15410157	50 µg	Hu	ChIP - ELISA - WB - IF
<a href="#">H3pan pAb</a>	C15410324	50 µg	Hu, BRE	ChIP - ELISA - WB - IF
<a href="#">H3pan mAb (clone B1B2)</a>	C15200011	50 µg	Hu, Mo, Ma, To, Po, Ar, BRE	ChIP - WB - IF
<a href="#">H3pan pAb</a>	C15310135	100 µl	Hu, Ze, BRE	ChIP - ELISA - WB
<a href="#">H4 pan pAb</a>	C15410156	50 µg	Hu, BRE	ChIP - ELISA - WB - IF
<a href="#">macroH2A.1/H2A.2 mAb</a>	C15210003	100 µg	Hu	ChIP-seq - ChIP - WB - IF
<a href="#">H2AX mAb</a>	C15210002	100 µg	Hu	ChIP - WB - IF
<a href="#">H2A.Z pAb</a>	C15410201	50 µg	Hu, Mo	ChIP-seq - ChIP - ELISA - WB - IF
<a href="#">H3.3 mAb</a>	C15210011	100 µg	Hu, BRE	ChIP-seq - ChIP - WB
<a href="#">H3K4un mAb</a>	C15200149	50 µg	Hu, BRE	ChIP - ELISA - WB - IF
<a href="#">H3K9un mAb</a>	C15200187	50 µg	Hu, BRE	ChIP - WB - IF

## Histone Acetylation



PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS
<a href="#">H2A.Zac pAb</a>	C15410202	50 µg	Hu, Mo	ChIP-seq - ChIP - ELISA - DB - WB - IF
<a href="#">H2A.Zac pAb</a>	C15410173	50 µg	Hu	ChIP-seq - ChIP - ELISA - DB - IF
<a href="#">H2A.ZK4ac mAb</a>	C15210009	100 µg	Hu	ChIP-seq - ChIP - WB - IF
<a href="#">H2A.ZK7ac mAb</a>	C15210012	100 µg	Hu	ChIP-seq - ChIP - WB - IF
<a href="#">H2AK5ac pAb</a>	C15410215	50 µg	Hu	ChIP-seq - ChIP - ELISA - DB - WB - IF
<a href="#">H2BK15ac pAb</a>	C15410220	50 µg	Hu	ChIP-seq - ChIP - ELISA - DB - WB - IF
<a href="#">H2BK12ac pAb</a>	C15410212	50 µg	Hu	ChIP-seq - ChIP - ELISA - DB - WB - IF
<a href="#">H2BK20ac mAb</a>	C15210010	100 µg	Hu	ChIP-seq - ChIP - WB - IF
<a href="#">H3K4ac pAb</a>	C15410322	50 µg	Hu, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
<a href="#">H3K9ac pAb</a>	C15410004	50 µg	Hu, Mo, Pi, Ze, Po, Pl, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
<a href="#">H3K9ac mAb</a>	C15200185	50 µg	Hu, BRE	ChIP - ELISA - WB - IF
<a href="#">H3K9acS10p pAb</a>	C15310102	100 µl	Hu, BRE	ChIP - ELISA - DB - WB - IF
<a href="#">H3K9/14ac pAb</a>	C15410200	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - PeA - WB - IF
<a href="#">H3K9/14ac pAb</a>	C15410005	50 µg	Hu, Mo, Ze, As, Ar, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
<a href="#">H3K14ac mAb</a>	C15210005	100 µg	Hu, BRE	ChIP-seq - ChIP - DB - WB - IF

pAb – polyclonal antibody mAb – monoclonal antibody

**APPLICATIONS:** ChIP – Chromatin immunoprecipitation ChIP-seq – ChIP followed by sequencing CUT&Tag – Cleavage under targets and tagmentation  
 ELISA - Enzyme-linked immunosorbent assay WB – Western blot DB – Dot blot IF – Immunofluorescence IS – Immunostaining PeA – Peptide array  
 PrA – Protein array IP - Immunoprecipitation FC - Flow cytometry

**REACTIVITY:** Al – Algae Ar – Arabidopsis As – Aspergillus Br – Brassica BRE – Broad range expected Co – Cow Dr – Drosophila Fu – Fungi  
 Hu – Human Ma – Maize Mo – Mouse Ne – Nematodes Pi – Pig Pl – Plasmodium Po – Poplar Ra – Rat Ri – Rice Sc – Schistosoma  
 St – Streptococcus To – Tomato TR – Trout Xe – Xenopus Ye – Yeast Ze – Zebrafish

PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS
H3K14ac pAb	C15410310	50 µg	Hu, BRE	ChIP - ELISA - DB - WB
H3K18ac pAb	C15410139	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
H3K23ac pAb	C15410344	50 µg	Hu, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
H3K27ac pAb	C15410196	50 µg	Hu, Mo, Ra, Ar, BRE	ChIP-seq - ChIP - ELISA - DB - PeA - WB - IF
H3K27ac pAb	C15410174	50 µg	Hu, Mo, Ra, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
H3K27ac mAb	C15200184	50 µg	Hu, BRE	ChIP - ELISA - WB - IF
H3K36ac pAb	C15410307	50 µg	Hu, BRE	ChIP-seq - ChIP - ELISA - DB - WB
H3K56ac pAb	C15410213	50 µg	Hu, BRE	ChIP-seq - ChIP - ELISA - DB
H3K79ac mAb	C15210006	100 µg	Hu, BRE	ChIP-seq - ChIP - WB - IF
H4K5ac pAb	C15410025	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
H4K5,8,12ac pAb	C15410021	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
H4K5,8,12,16ac pAb	C15410024	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - IF
H4K8ac pAb	C15410103	50 µg	Hu, Mo, BRE	ChIP - ELISA - DB - WB - IF
H4K8ac mAb	C15200217	50 µg	Hu, BRE	ChIP - DB - WB
H4K12ac pAb	C15410331	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
H4K12ac mAb	C15200218	50 µg	Hu, BRE	ChIP - WB
H4K16ac mAb	C15200219	50 µg	Hu, BRE	ChIP - DB - WB
H4K20ac mAb	C15210008	100 µg	Hu, BRE	ChIP-seq - ChIP - WB - IF

## Histone methylation

SAMPLE SIZE  
AVAILABLE

PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS
H3K4me1 pAb	C15410194	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - PeA - WB - IF
H3K4me1 pAb	C15410037	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
H3K4me1 mAb	C15200150	50 µg	Hu, BRE	ChIP - ELISA - WB - IF
H3K4me2 pAb	C15410035	50 µg	Hu, Ar, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
H3K4me2 mAb	C15200151	50 µg	Hu, As, BRE	ChIP - ELISA - WB - IF
H3K4me3 pAb	C15410003	50 µg	Hu, Mo, Ar, Ze, Tr, Ri, To, Ma, Po, BRE	ChIP-seq - ChIP - CUT&Tag ELISA - DB - PeA - WB - IF
H3K4me3 pAb	C15410030	50 µg	Hu, Mo, Ar, BRE	ChIP-seq - ChIP - DB - WB - IF
H3K4me3 mAb	C15200152	50 µg	Hu, Mo, Ne, Ar, BRE	ChIP-seq - ChIP - ELISA - WB - IF
H3K4me3 pAb	C15310003	100 µl	Hu, Ne, BRE	ChIP - ELISA - DB - WB
H3K9me1 pAb	C15410065	50 µg	Hu, BRE	ChIP - ELISA - DB - WB
H3K9me2 pAb	C15410060	50 µg	Hu, Mo, Ar, Ri, To, Br, BRE	ChIP - ELISA - DB - WB - IF
H3K9me2 mAb	C15200154	50 µg	Hu, Ze, Ar, ,BRE	ChIP - ELISA - WB - IF
H3K9me3 pAb	C15410193	50 µg	Hu, Mo, Ye, BRE	ChIP-seq - ChIP - CUT&Tag- ELISA - DB - PeA - WB - IF
H3K9me3 pAb	C15410056	50 µg	Hu, Mo, Ze, Tr, Ar, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
H3K9me3 rAb	C15500003	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - IF



Click on the antibody's name to visit the product page

PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS
<a href="#">H3K9me3 mAb</a>	C15200146	50 µg	Hu, Mo, Fu, BRE	ChIP - DB - WB - IF
<a href="#">H3K9me3 mAb</a>	C15200153	50 µg	Hu, Fu, BRE	ChIP - ELISA - WB - IF
<a href="#">H3K9me3S10p pAb</a>	C15310128	100 µl	Hu, BRE	ChIP - ELISA - DB - WB - IF
<a href="#">H3K23me2 mAb</a>	C15210007	100 µg	Hu, BRE	ChIP-seq - ChIP - WB
<a href="#">H3K27me1 pAb</a>	C15410045	50 µg	Hu, Ar, BRE	ChIP - ELISA - DB - WB - IF
<a href="#">H3K27me2 pAb</a>	C15410046	50 µg	Hu, Ze, BRE	ChIP - ELISA - DB - WB - IF
<a href="#">H3K27me3 pAb</a>	C15410195	50 µg	Hu, Mo, Dr, Ne, Ar, Ma, To, Po, Al, BRE	ChIP-seq - ChIP - CUT&Tag - ELISA - DB - WB - IF
<a href="#">H3K27me3 pAb</a>	C15410069	50 µg	Hu, Mo, Ra, Pi, Co, Ze, Dr, Sc, Ar, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
<a href="#">H3K27me3 mAb</a>	C15200181	50 µg	Hu, BRE	ChIP - ELISA - WB - IF
<a href="#">H3K36me1 pAb</a>	C15410089	50 µg	Hu, BRE	ChIP - ELISA - DB - WB - IF
<a href="#">H3K36me2 pAb</a>	C15310127	100 µl	Hu, Mo, Ye, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
<a href="#">H3K36me2 mAb</a>	C15200182	50 µg	Hu, BRE	ChIP - ELISA - DB - WB
<a href="#">H3K36me3 pAb</a>	C15410192	50 µg	Hu, Mo, Ar, Ri, BRE	ChIP-seq - ChIP - ELISA - DB - PeA - WB - IF
<a href="#">H3K36me3 pAb</a>	C15410058	50 µg	Hu, Ze, Dr, BRE	ChIP-seq - ChIP - ELISA - DB - WB
<a href="#">H3K36me3 mAb</a>	C15200183	50 µg	Hu, Ra, BRE	ChIP - ELISA - DB - WB - IF
<a href="#">H3K79me1 pAb</a>	C15410082	50 µg	Hu, Ye, BRE	ChIP-seq - ChIP - ELISA - DB - WB
<a href="#">H3K79me2 pAb</a>	C15410051	50 µg	Hu, Yet, BRE	ChIP-seq - ChIP - ELISA - DB - WB
<a href="#">H3K79me2 pAb</a>	C15310051	100 µl	Hu, BRE	ChIP - ELISA - DB - WB
<a href="#">H3K79me3 pAb</a>	C15410068	50 µg	Hu, Mo, Yet, BRE	ChIP-seq - ChIP - ELISA - DB - WB
<a href="#">H3R17me2(asym)K18ac pAb</a>	C15410171	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF
<a href="#">H4K20me1 mAb</a>	C15200147	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - DB - WB - IF
<a href="#">H4K20me1 pAb</a>	C15410034	50 µg	Hu, BRE	ChIP - ELISA - DB - WB - IF
<a href="#">H4K20me2 mAb</a>	C15200220	50 µg	Hu, BRE	ChIP - DB - WB - IF
<a href="#">H4K20me3 pAb</a>	C15410207	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - PeA - WB - IF
<a href="#">H4K20me3 pAb</a>	C15410057	50 µg	Hu, Mo, BRE	ChIP-seq - ChIP - ELISA - DB - WB - IF

## Histone phosphorylation

SAMPLE SIZE  
AVAILABLE

PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS
<a href="#">H3S10p pAb</a>	C15410116	50 µg	Hu, BRE	ChIP - ELISA - DB - WB - IF
<a href="#">H3T3p mAb</a>	C15210001	100 µg	Hu, BRE	ChIP - WB - IF
<a href="#">H3T11p pAb</a>	C15410309	50 µg	Hu, BRE	ChIP - ELISA - DB - WB - IF
<a href="#">H4S1p mAb</a>	C15210004	100 µg	Hu, BRE	ChIP - WB - IF

# Histone ubiquitination

SAMPLE SIZE  
AVAILABLE

PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS
<a href="#">H2AK119ub pAb</a>	C15410002	50 µg	Hu	ChIP - ELISA - DB - WB
<a href="#">yeast H2BK123ub mAb</a>	C15200232	50 µg	Hu, Ye	ChIP-seq - ChIP - WB

# Negative control IgG

PRODUCT NAME	CAT. NO.	FORMAT
<a href="#">Rabbit IgG</a>	C15410206	250 µg
<a href="#">Mouse IgG</a>	C15400001	100 µg
<a href="#">Rat IgG</a>	C15420001	25 µg

# Non-histone proteins

## Transcription

SAMPLE SIZE  
AVAILABLE

PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS	siRNA VALIDATED
<a href="#">AF9 pAb</a>	C15310266	100 µl	Hu	ChIP-seq - ChIP - ELISA	
<a href="#">AML1-ETO pAb</a>	C15310197	100 µl	Hu	ChIP-seq - ChIP - ELISA - WB	
<a href="#">CBFb pAb</a>	C15310002	100 µl	Hu	ChIP-seq - ChIP - ELISA	
<a href="#">CHD7 pAb</a>	C15410340	100 µl	Hu	ChIP-seq - ChIP - WB - IP	
<a href="#">CTCF pAb</a>	C15410210	50 µg	Hu, Mo	ChIP-seq - ChIP - ELISA - WB - IF	✓
<a href="#">EHF pAb</a>	C15410363	50 µg	Hu	ChIP-seq - ChIP - ELISA	
<a href="#">ER alpha mAb</a>	C15100066	100 µl	Hu, Mo	ChIP-seq - ChIP - ELISA - WB - IHC - IP	
<a href="#">ER alpha mAb</a>	C15200009	50 µg	Hu	ChIP - WB - IHC	
<a href="#">ETO pAb</a>	C15310001	100 µl	Hu	ChIP-seq - ChIP - ELISA	
<a href="#">FOXA2 pAb</a>	C15410343	50 µg	Hu	ChIP-seq - ChIP - ELISA - WB	
<a href="#">GR mAb</a>	C15200010	50 µg	Hu, Ra	ChIP-seq - ChIP - ELISA - WB - IHC - FC - IP	
<a href="#">NRF1 mAb</a>	C15200013	50 µg	Hu	ChIP-seq - ChIP - WB	
<a href="#">OCT4 pAb</a>	C15410305	50 µg	Hu	ChIP-seq - ChIP - ELISA - WB	
<a href="#">p53 pAb</a>	C15410083	50 µg	Hu	ChIP-seq - ChIP - ELISA - WB	
<a href="#">Pol II mAb</a>	C15100055	100 µl	Hu, Mo, Xe	ChIP-seq - ChIP - WB	✓
<a href="#">Pol II mAb</a>	C15200004	50 µg	Hu	ChIP-seq - ChIP - ELISA - WB - IF	✓
<a href="#">Pol II S2p mAb</a>	C15200005	50 µg	Hu, Ze	ChIP-seq - ChIP - ELISA - WB - IF	✓
<a href="#">Pol II S5p mAb</a>	C15200007	50 µg	Hu	ChIP-seq - ChIP - ELISA - WB - IF	✓
<a href="#">PPARG pAb</a>	C15410367	50 µg	Hu	ChIP-seq - ChIP - ELISA	



Click on the antibody's name to visit the product page

PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS	siRNA VALIDATED
<a href="#">RARA pAb</a>	C15310155	100 µl	Hu	ChIP-seq - ChIP - ELISA - WB	
<a href="#">TAL1 mAb</a>	C15200012	50 µg	Hu	ChIP-seq - ChIP - IF	
<a href="#">TBP mAb</a>	C15200002	100 µg	Hu, Mo	ChIP-seq - ChIP - WB	✓
<a href="#">YY1 pAb</a>	C15410345	50 µg	Hu	ChIP-seq - ChIP - ELISA - WB	
<a href="#">ZMYM3 mAb</a>	C15200016	50 µg	Hu	ChIP-seq - ChIP - PrA - IF	

## Chromatin associated protein

SAMPLE SIZE  
AVAILABLE

PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS	siRNA VALIDATED
<a href="#">BRD4 pAb</a>	C15410337	50 µg	Hu	ChIP-seq - ChIP - ELISA - WB	
<a href="#">CHD1 pAb</a>	C15410334	100 µl	Hu	ChIP-seq - ChIP - WB - IP	
<a href="#">CXXC1 pAb</a>	C15410315	50 µg	Hu	ChIP-seq - ChIP - WB - IF	
<a href="#">MeCP2 pAb</a>	C15410052	50 µg	Hu	ChIP - ELISA - WB	✓
<a href="#">NCOR1 pAb</a>	C15410341	100 µl	Hu, Mo	ChIP-seq - ChIP - WB - IP	
<a href="#">p300 mAb</a>	C15200211	50 µg	Hu	ChIP-seq - ChIP	
<a href="#">RBBP5 pAb</a>	C15410342	100 µl	Hu	ChIP-seq - ChIP - WB - IP - IHC	
<a href="#">TIP-5 pAb</a>	C15310090	100 µl	Hu, Mo	ChIP - WB	

# Histone modifying enzymes

SAMPLE SIZE  
AVAILABLE

PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS	siRNA VALIDATED
<a href="#">CBX2 pAb</a>	C15410339	100 µl	Hu	ChIP-seq - ChIP - WB - IP	
<a href="#">CBX8 pAb</a>	C15410333	100 µl	Hu	ChIP-seq - ChIP - WB - IP	
<a href="#">EZH2 pAb</a>	C15410039	50 µg	Hu, Mo	ChIP-seq - ChIP - WB - IF	✓
<a href="#">EZH2 mAb</a>	C15200180	50 µg	Hu, Mo	ChIP	
<a href="#">HDAC1 pAb</a>	C15410325	50 µg	Hu, Mo	ChIP-seq - ChIP - ELISA - WB - IF - PrA	✓
<a href="#">HDAC2 mAb</a>	C15200201	50 µg	Hu	ChIP-seq - ChIP - WB - IF	✓
<a href="#">HDAC3 pAb</a>	C15410364	50 µg	Hu	ChIP-seq - ChIP - ELISA	
<a href="#">HDAC6 pAb</a>	C15410362	50 µg	Hu	ChIP-seq - ChIP - ELISA	
<a href="#">JARID1C pAb</a>	C15410338	100 µl	Hu, Mo	ChIP-seq - ChIP - WB - IP	
<a href="#">JMJD2a pAb</a>	C15410332	100 µl	Hu, Mo	ChIP-seq - ChIP - WB - IP	
<a href="#">KAT2B pAb</a>	C15410335	100 µl	Hu	ChIP-seq - ChIP - WB	
<a href="#">LSD1 pAb</a>	C15410067	50 µg	Hu	ChIP-seq - ChIP - ELISA - WB - IF	✓
<a href="#">MLL1 pAb</a>	C15310264	100 µl	Hu	ChIP-seq - ChIP - ELISA	
<a href="#">PHF8 pAb</a>	C15410336	100 µl	Hu	ChIP - WB - IP	
<a href="#">SAP30 pAb</a>	C15410036	50 µg	Hu	ChIP-seq - ChIP - WB	
<a href="#">WDR5 pAb</a>	C15410027	50 µg	Hu	ChIP-seq - ChIP - WB - IF	

## Other

SAMPLE SIZE  
AVAILABLE

PRODUCT NAME	CAT. NO.	FORMAT	REACTIVITY	APPLICATIONS
<a href="#">CRISPR/Cas9 pAb</a>	C15310258	100 µl	St	ChIP-seq - ChIP - WB - IF - IP
<a href="#">CRISPR/Cas9 mAb</a>	C15200229	50 µg	St	ChIP - WB - IF - IP
<a href="#">CRISPR/Cas9 C-terminal mAb</a>	C15200231	50 µg	St	ChIP - WB - IF
<a href="#">HA tag mAb</a>	C15200190	50 µg	NA	ChIP - IP - WB - IS
<a href="#">MYH11 pAb</a>	C15310254	100 µl	Hu	ChIP-seq - ChIP - ELISA - WB
<a href="#">Ty1 mAb</a>	C15200054	50 µg	NA	ChIP - WB

## Negative control IgG

PRODUCT NAME	CAT. NO.	FORMAT
<a href="#">Rabbit IgG</a>	C15410206	250 µg
<a href="#">Mouse IgG</a>	C15400001	100 µg
<a href="#">Rat IgG</a>	C15420001	25 µg

Shop online in our EpiStore at  
**[www.diagenode.com](http://www.diagenode.com)**