



HOLOGIC®
Diagenode

**Plant Epigenetics
Antibodies**

About us

With over 20 years of expertise in epigenomics, Hologic Diagenode is your trusted supplier of **validated, high-quality antibodies**.

Why Choose Us?



High Quality Antibodies

Antibodies validated for epigenetic applications like ChIP, ChIP-seq, and CUT&Tag, ensuring high specificity and reproducibility.



Performance Guarantee

We conduct strict in-house validations and lot-to-lot comparisons, guaranteeing the performance of antibodies for tested applications.



Transparency and Accessibility

We provide batch-specific validation data on our website, ensuring transparency and accessibility for researchers.

Antibody	Cat. No.	Format	Applications tested
5-methylcytosine (5-mC)	C15200006	100 µg	MeDIP - IF
H3pan 1B1B2	C15200011	50 µg	ChIP - WB - IF
H3K4me3	C15200152	50 µg	ChIP/ChIP-seq - CUT&Tag - ELISA - WB - IF
H3K9me2	C15200154	50 µg	ChIP - ELISA - WB - IF
H3K4me3	C15410003	50 µg	ChIP/ChIP-seq - ELISA - Dot Blot - WB - IF
H3K9ac	C15410004	50 µg	ChIP/ChIP-seq - ELISA - Dot Blot - WB - IF
H3K9/14ac	C15410005	50 µg	ChIP/ChIP-seq - ELISA - Dot Blot - WB - IF
H3K4me3	C15410030	50 µg	ChIP/ChIP-seq - Dot Blot - WB - IF
H3K4me2	C15410035	50 µg	ChIP/ChIP-seq - CUT&Tag - ELISA - Dot Blot - WB
H3K27me1	C15410045	50 µg	ChIP - ELISA - Dot Blot - WB - IF
H3K9me3	C15410056	50 µg	ChIP/ChIP-seq - ELISA - Dot Blot - WB
H3K9me2	C15410060	50 µg	ChIP - ELISA - Dot blot - WB - IF
H3K27me3	C15410069	50 µg	ChIP/ChIP-seq - ELISA - Dot blot - WB - IF
H3K36me3	C15410192	50 µg	ChIP/ChIP-seq - CUT&Tag - ELISA - Dot blot - peptide array - WB - IF
H3K27me3	C15410195	50 µg	ChIP/ChIP-seq - CUT&Tag - ELISA - Dot blot - WB - IF
H3K27ac	C15410196	50 µg	ChIP/ChIP-seq - CUT&Tag - ELISA - Dot blot - WB - IF
H3R2me1	C15410276	50 µg	WB
H3T3pK4me1	C15410277	50 µg	Dot blot - WB - IF - IHC
H3T3pK4me2	C15410278	50 µg	ChIP - Dot blot - WB - IF - IHC
H3T3pK4ac	C15410279	50 µg	ChIP - Dot blot - WB - IF - IHC
H3K4me1T6p	C15410280	50 µg	ChIP - Dot blot - WB - IF - IHC
H3K4me3T6p	C15410281	50 µg	ChIP - Dot blot - WB - IF - IHC
H3T6p	C15410282	50 µg	ChIP - Dot blot - WB - IF - IHC
H3T6pK9me1	C15410283	50 µg	ChIP - Dot blot - WB - IF - IHC
H3T6pK9me2	C15410284	50 µg	ChIP - WB - IF - IHC
H3T6pK9me3	C15410285	50 µg	ChIP - Dot blot - WB - IF - IHC
H3R8me2(asym)	C15410286	50 µg	ChIP - WB - IF - IHC
H3R8me2(sym)	C15410287	50 µg	ChIP - WB
H3S10pT11p	C15410288	50 µg	ChIP - WB - IF - IHC
H3R17me2(asym)	C15410289	50 µg	ChIP - WB - IF
H3K18me1	C15410290	50 µg	ChIP - Dot blot - WB - IF - IHC
H3K18me2	C15410291	50 µg	ChIP - Dot blot - WB - IF - IHC
H3K27me3S28p	C15410293	50 µg	ChIP - Dot blot - WB - IF - IHC
H3R2me2K4me2	C15410294	50 µg	ChIP - Dot blot - WB - IF - IHC
H3K37me1	C15410295	50 µg	Dot blot - WB
H3K56me1	C15410296	50 µg	ChIP - Dot blot - WB - IF
H3K56me3	C15410297	50 µg	ChIP - Dot blot - WB - IF
H4K16ac	C15410300	50 µg	Dot blot - WB - IF

Contact Us

EU // Liège Science Park / Rue Bois Saint-Jean 3, 4102 Seraing / BE / Phone : +32 4 364 20 50 / Fax : +32 4 364 20 51

US // 400 Morris Avenue / Suite 101 Denville / NJ 07834 / USA / Phone : +1 862 209 4680 / Fax : +1 862 209 4681

www.diagencode.com