



HOLOGIC[®] Diagenode

Plant Epigenetics Kits & Reagents

About us

With over 20 years in epigenomics, Hologic Diagenode is your trusted partner for **high-quality reagents** and **kits**. We are committed to excellence, providing reliable tools to advance your research with confidence.



Chromatin

Your all-in-one solution for every step of the protocol, validated and proven effective across diverse plant species.

Universal Plant ChIP-seq Kit

Cat. No. C01010152



Streamline Library Preparation with our Optimized Kit.

MicroPlex Library Preparation Kit v3

Cat. No. C05010001



Optimizing the Assessment of Open Chromatin Regions.

ATAC-seq Kit

Cat. No. C01080001





DNA Methylation

Genome-wide studies at single nucleotide resolution with modified RRBS

Gene-specific or genome-wide studies at 100–500 bp resolution with MeDIP

Premium RRBS kit V2

Cat. No. C02030036



MagMeDIP Kit

Cat. No. C02010021



Small RNAs

Enhance your transcriptomic workflow by integrating the D-Plex Small RNA kit, facilitating the discovery of novel small non-coding RNAs and optimizing ribosomal profiling.

D-Plex Small RNA-seq Kit for Illumina

Cat. No. C05030001



Tagmentation-Based Approaches

Leverage our Tagmentase (Tn5 transposase) for all your tagmentation-based applications.

Tagmentase (Tn5 transposase) – loaded

Cat. No. C01070012



Tagmentase (Tn5 transposase) - unloaded

Cat. No. C01070010



Tagmentation Buffer (2x)

Cat. No. C01019043



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.diagenode.com