

DNA METHYLATION

GENE-SPECIFIC

Immunoprecipitation

- All species
- Methylation status of specific region

MeDIP-qPCR

hMeDIP-qPCR

MethylCap-qPCR
(MBD capture)

GENOME-WIDE

Immunoprecipitation

- All species
- All methylated regions

MeDIP-seq

MethylCap-seq
(MBD capture)

REDUCED GENOME

Bisulfite conversion

- All vertebrates
- 4 M CpGs (>10x)

RRBS

Reduced Representation
Bisulfite Sequencing

100-500 bp resolution

Single-base resolution

PRODUCTS

KITS

Methylated DNA immunoprecipitation	Cat. No.	Format
MagMeDIP kit	C02010020	10 rxns
	C02010021	48 rxns
DNA methylation control package	C02040012	40 rxns
DNA methylation control package V2	C02040019	48 rxns
MethylCap kit	C02020010	48 rxns
MethylCap protein	C02020012	100 µg
Hydroxymethylated DNA immunoprecipitation		
hMeDIP kit	C02010031	16 rxns
Bisulfite conversion		
Premium RRBS kit V2	C02030036	24 rxns
	C02030037	96 rxns
Premium Methyl UDI-UMI Adapters Module - Set A	C02030040	24 rxns
Premium Methyl UDI-UMI Adapters Module - Set B	C02030041	24 rxns
Premium Bisulfite kit	C02030030	50 rxns

Antibodies

5-methylcytosine (5-mC) (clone 33D3)	C15200081	DNMT1	C15310111
5-hydroxymethylcytosine (5-hmC) (mouse)	C15200200	DNMT3A	C15410084
3-methylcytosine (3-mC)	C15410209	DNMT3B	C15410218
5-Carboxylcytosine (5-caC)	C15410204	TDG	C15200227
5-formylcytosine (5-fC)	C15310200	TET2	C15200179