

DNA hydroxymethylation control package

Cat. No. **C02040018**

Format: 48 rxns

Product content

Component	Quantity	Concentration	Storage
Hydroxymethylated spike-in control*	48 µl	0.001 ng/µl	-20°C/-4°F
Primer pair for hydroxymethylated spike-in ctrl**	144 µl	5 µM each	-20°C/-4°F

* Alternatively named hmeDNA control in the hMeDIP kit (Cat. No. C02010031)

** Alternatively named hmeDNA primer pairs in the hMeDIP kit (Cat. No. C02010031)

Product description

The DNA hydroxymethylation control package includes one hydroxymethylated spike-in control and its corresponding qPCR primer set that can be added to the DNA sample of interest for any hydroxymethylation profiling experiment.

The hydroxymethylated spike-in control is produced from a genomic sequence from *Arabidopsis thaliana* and may therefore interfere with plant samples. However, it does not show significant homology with other samples species (e.g. human, mouse or rat).

Applications

For capture of hydroxymethylated DNA (e.g. hMeDIP - Hydroxymethylated DNA Immunoprecipitation)

The spike-in control must be added to the DNA sample of interest before the capture. The package also includes a primer pair targeting the hydroxymethylated spike-in control, allowing to check the success of the capture by qPCR. The size of the expected amplicon is 88 bp.

Storage conditions: Store at -20°C/-4°F. For long storage, store at -80°C /-112°F. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: December 16, 2020