
5-Hydroxymethylcytosine Antibody [HMC/4D9]

(Cat. No. A-1018)

Background

5-hydroxymethylcytosine (5-hmC), as a sixth DNA base with functions in transcription regulation, has been detected to be abundant in human and mouse brain and embryonic stem (ES) cells. In mammals, it can be generated by oxidation of 5-mC, a reaction mediated by the ten-eleven translocation (TET) family of 5mC-hydroxylases. 5-hmC was demonstrated to play an important and different role than 5-mC in the regulation of DNA methylation, chromatin remodeling, and gene expression.

Concentration

1 mg/ml

Description

Mouse monoclonal antibody to 5-Hydroxymethylcytosine (5-hmC), clone HMC/4D9

Specificity

Modified base 5-hydroxymethylcytosine (5-hmC), a broad range of species

Isotype

IgG1

Formulation

10 mM phosphate buffer, 150 mM NaCl, pH 7.4.

Storage

4°C. For long-term storage, aliquot and store at -20°C. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance. Stable for 6 months from the date of shipment.

Purity

Protein A purified

Handling Recommendations

For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Alternate Names

5-hydroxymethylcytidine, anti-5-hydroxymethylcytidine, anti-5-hydroxymethylcytosine, 5hmC, 5hmeC, 5-hmeC

Application

Immunofluorescence: 1:200 – 1: 500

Immunohistochemistry: 1:200 – 1: 500

ELISA: 1:1000 – 1:2000

Dot Blot: 1:2000

hMeDIP: 0.5-1 µg/10⁶ Cells

Ordering Information

Products	Size	Cat. No.
5-Hydroxymethylcytosine Antibody [HMC/4D9]	10 µg	A-1018-010
	50 µg	A-1018-050
	100 µg	A-1018-100

This product is for research purposes only. Not intended for use in diagnostic procedures.