

PRODUCT INFORMATION

Anti-5-hydroxy Methylcytosine Rabbit Monoclonal Antibody

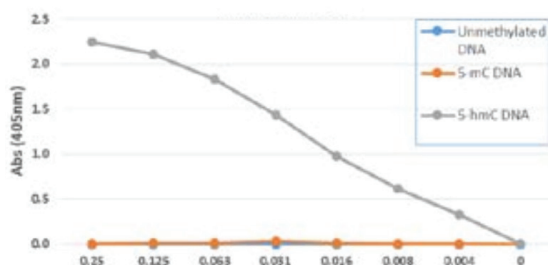
(Clone RM236)

Item No. 20723

Overview and Properties

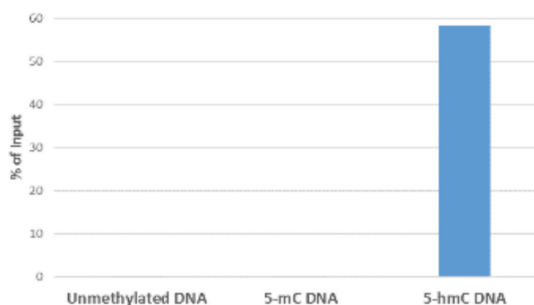
Contents:	This vial contains 50 µg of protein A affinity-purified antibody from an animal origin free culture supernatant.
Synonym:	5-hmC
Immunogen:	BSA-conjugated 5-hydroxy methylcytosine
Cross Reactivity:	(+) 5-hydroxy methylcytosine in both single-stranded and double-stranded DNA; (-) Non-methylated cytosine, methylcytosine in DNA
Species Reactivity:	(+) All species
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	As supplied, 1 year from the QC date provided on the Certificate of Analysis, when stored properly
Storage Buffer:	50% glycerol/PBS with 1% BSA and 0.09% sodium azide
Clone:	RM236
Host:	Rabbit
Isotype:	IgG
Applications:	Dot blot (DB), ELISA, Hydroxymethylated DNA immunoprecipitation (hMeDIP), Immunocytochemistry (ICC), and Immunohistochemistry (IHC); the recommended starting dilution for DB is 0.2 - 1 µg/ml, 0.1 - 1 µg/ml for ELISA and IHC, 0.2 - 2 µg/ml for hMeDIP, and 0.5 - 2 µg/ml for ICC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Anti-5-hydroxy Methylcytosine Rabbit Monoclonal Antibody (Clone RM236) (µg/ml)

ELISA of single-stranded DNA using Anti-5-hydroxy Methylcytosine Rabbit Monoclonal Antibody (Clone RM236). The plate was coated with streptavidin and then biotinylated single-stranded unmethylated DNA, 5-methylcytosine (5-mC) DNA, and 5-hydroxy methylcytosine (5-hmC) DNA. A serial dilution of RM236 was used as the primary antibody, and an alkaline phosphatase conjugated anti-rabbit IgG as the secondary antibody.



hMeDIP was performed using anti-5-hmC antibody (RM236) at a 10:1 DNA:Ab ratio. 1 ng of unmethylated, 5-methylcytosine (5-mC) or 5-hydroxy methylcytosine (5-hmC) DNA standard (897 bp) was spiked in 1 µg of genomic DNA isolated from HeLa cells as the control. Realtime PCR was then performed to determine the capture of DNA standard as in % of input.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

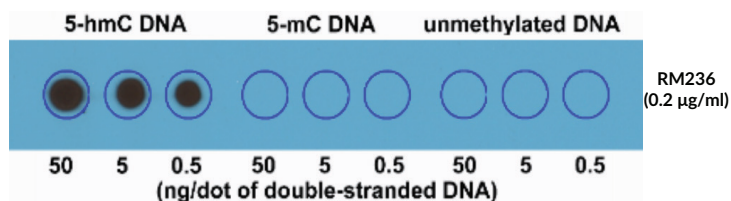
SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

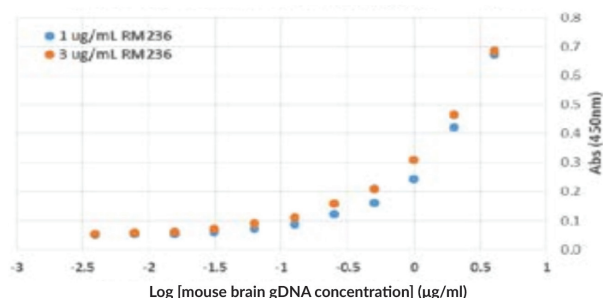
Copyright Cayman Chemical Company, 11/11/2020

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

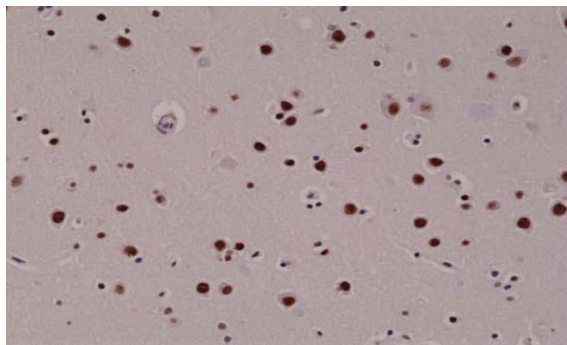
PRODUCT INFORMATION



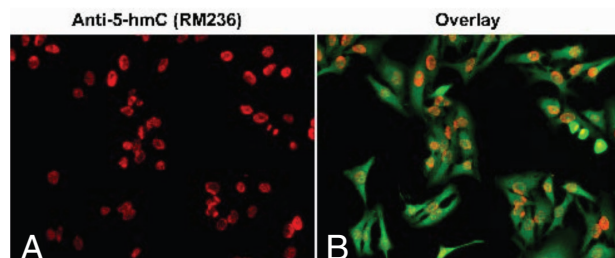
Dot blot of double-stranded DNA using Anti-5-hydroxy Methylcytosine Rabbit Monoclonal Antibody (Clone RM236). The membrane was pre-spotted with 50, 5, and 0.5 ng/dot of double stranded 5-hydroxy methylcytosine (5-hmC) DNA, 5-methylcytosine (5-mC) DNA, and unmethylated DNA.



Direct ELISA of mouse brain genomic DNA using anti-5-hmC antibody (RM236). The plate was directly coated with different concentrations of genomic DNA isolated from mouse brain tissue. 1 µg/ml or 3 µg/ml of RM236 was used as the primary antibody, and a HRP conjugated anti-rabbit IgG as the secondary antibody.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human brain tissue sections, using Anti-5-hydroxy Methylcytosine Rabbit Monoclonal Antibody (Clone RM236).



Panel A: Immunocytochemical labeling of HeLa cells using 0.5 µg/ml Anti-5-hydroxy Methylcytosine Rabbit Monoclonal Antibody (Clone RM236) (red). Panel B: Actin filaments were labeled with fluorescein phalloidin (green). HeLa cells were fixed with 4% paraformaldehyde and permeabilized with methanol (-20°C) before treatment with 2 N HCl for 30 minutes at 37°C to denature the DNA.

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

5-Hydroxymethylcytosine (5-hmC) is an epigenetic modification formed from the oxidation of 5-methylcytosine by Tet dioxygenases.¹ 5-hmC is primarily a stable DNA modification, but it can be oxidized by Tet enzymes and its products further modified to generate nonmethylated cytosine, indicating a role as an intermediate in DNA demethylation as well.²⁻⁴ It is associated with actively transcribed genes and recognized by a variety of proteins, including proteins involved in DNA repair and DNA metabolic processes.^{2,3} 5-hmC has been found in all mammalian tissues but levels are higher in the brain relative to other tissues.⁵ The percentage of genomic 5-hmC in mouse cerebellum increases during postnatal development until adulthood, and genes acquiring intragenic 5-hmC are enriched in pathways associated with age-related neurodegenerative disease pathways in adult mice.⁶ In contrast, 5-hmC levels are reduced by up to 8-fold in cancer tissues.^{1,7} Cayman's 5-Hydroxymethylcytosine Polyclonal Antibody can be used for dot blot, ELISA, hydroxymethylated DNA immunoprecipitation (hMeDIP), immunocytochemistry (ICC), and immunohistochemistry (IHC) applications.

Reference

1. Tahiliani, M., Koh, K.P., Shen, Y., *et al.* Conversion of 5-methylcytosine to 5-hydroxymethylcytosine in mammalian DNA by MLL partner TET1. *Science* **324**(5929), 930-935 (2009).
2. Bachman, M., Uribe-Lewis, S., Yang, X., *et al.* 5-Hydroxymethylcytosine is a predominantly stable DNA modification. *Nat. Chem.* **6**(12), 1049-1055 (2014).
3. Spruijt, C.G., Gnerlich, F., Smits, A.H., *et al.* Dynamic readers for 5-(hydroxy)methylcytosine and its oxidized derivatives. *Cell* **152**(5), 1146-1159 (2013).
4. Breiling, A. and Lyko, F. Epigenetic regulatory functions of DNA modifications: 5-methylcytosine and beyond. *Epigenetics Chromatin* **8**, 24 (2015).
5. Globisch, D., Münzel, M., Müller, M., *et al.* Tissue distribution of 5-hydroxymethylcytosine and search for active demethylation intermediates. *PLoS One* **5**(12), e15367 (2010).
6. Song, C.X., Szulwach, K.E., Fu, Y., *et al.* Selective chemical labeling reveals the genome-wide distribution of 5-hydroxymethylcytosine. *Nat. Biotechnol.* **29**(1), 68-72 (2011).
7. Wang, Z., Du, M., Yuan, Q., *et al.* Epigenomic analysis of 5-hydroxymethylcytosine (5hmC) reveals novel DNA methylation markers for lung cancers. *Neoplasia* **22**(3), 154-161 (2020).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM