

PRODUCT INFORMATION



5-Methylcytosine DNA Matched Antibody Pair (Clones RM231/RM476)

Item No. 42873

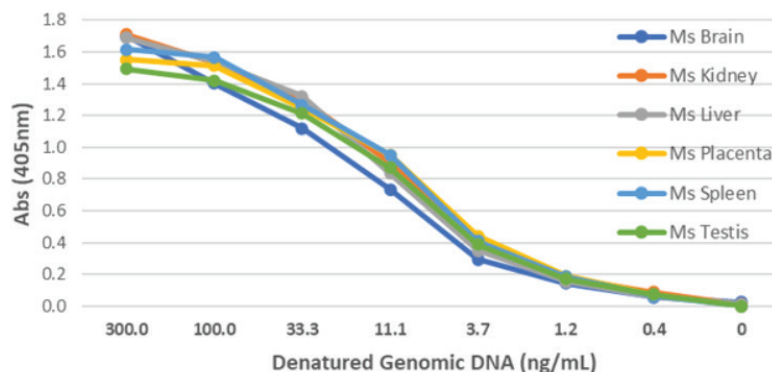
Overview and Properties

Antibody	Item Name	Storage Buffer	Clone	Applications*
Capture	5-Methylcytosine Rabbit Monoclonal Antibody (Clone RM231) (Item No. 20722)	50 µg at 1.0 mg/ml in PBS with 50% glycerol and 0.09% sodium azide	RM231	Sandwich ELISAs: 50-200 ng/well
Detection	Adenine and Guanine Rabbit Monoclonal Antibody - Biotinylated (Clone RM476)	25 µg at 1.0 mg/ml in PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide	RM476	Sandwich ELISAs: 0.1-0.5 µg/ml

*Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Contents: Contains protein A-affinity purified monoclonal antibodies.
Synonyms: 5-mC, m⁵C, NSC 137776
Immunogen: BSA-conjugated 5-mC
Cross Reactivity: (+) 5-mC; (-) Non-methylated cytosine, 5-hmC, 5-caC, 5-fC
Form: Liquid
Storage: -20°C (as supplied)
Stability: ≥1 year
Host: Rabbit

Image



Detection of 5-methylcytosine in genomic DNA from different mouse tissues. Results are shown for a sandwich ELISA using 5-Methylcytosine Rabbit Monoclonal Antibody (RM231) (Item No. 20722) as the capture antibody (50 ng/well at 1 µg/ml) and Adenine-Guanine Rabbit Monoclonal Antibody - Biotinylated (Clone RM476) as the detection antibody (0.2 µg/ml), followed by AP-conjugated streptavidin.

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, 07/01/2026

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-Methylcytosine (5-mC) is an epigenetic modification formed from methylation of cytosine by DNA methyltransferases (DNMTs).¹ 5-mC is an abundant DNA modification that is enriched at CpG dinucleotides. It can be successively oxidized by ten-eleven translocation (TET) dioxygenases to form 5-hydroxymethylcytosine (5-hmC), 5-formylcytosine (5-fC), and 5-carboxylcytosine (5-caC). 5-mC is associated with reduced gene transcription of DNA but is also found in the tRNA, mRNA, and rRNA from bacteria, archaea, and eukaryotes.^{1,2} Levels of 5-mC on CpG islands in the genes encoding p16, thrombospondin-1 (TSP-1), and methylated in tumor 1 (MINT1) are increased in tumors compared with adjacent normal tissue in patients with renal cell carcinoma.³ Cayman's 5-Methylcytosine DNA Matched Antibody Pair (Clones RM231/RM476) contains 5-Methylcytosine Rabbit Monoclonal Antibody (Clone RM231) (Item No. 20722) as a capture antibody and Adenine-Guanine Rabbit Monoclonal Antibody - Biotinylated (Clone RM476) as a detection antibody. This antibody pair can be used for sandwich ELISAs or other sandwich assays for the specific detection of 5-mC in ssDNA.

References

1. Breiling, A. and Lyko, F. Epigenetic regulatory functions of DNA modifications: 5-Methylcytosine and beyond. *Epigenetics Chromatin* **8**, 24 (2015).
2. Edelheit, S., Schwartz, S., Mumbach, M.R., *et al.* Transcriptome-wide mapping of 5-methylcytidine RNA modifications in bacteria, archaea, and yeast reveals m⁵C within archaeal mRNAs. *PLoS Genetics* **9**(6), e1003602 (2013).
3. Arai, E., Kanai, Y., Ushijima, S., *et al.* Regional DNA hypermethylation and DNA methyltransferase (DNMT) 1 protein overexpression in both renal tumors and corresponding nontumorous renal tissues. *Int. J. Cancer* **119**(2), 288-296 (2006).

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