

5-Methylcytosine (5-Methylcytidine) Monoclonal Antibody [33D3]

(Catalog No. A-1014)

Background:	DNA methylation is the major modification of eukaryotic genomes and plays an essential role in mammalian development. DNA methylation of promoter regions leads to inactivation of gene function. Also, DNA methylation varies according to tissue type, and Region-specific DNA hypermethylation and global DNA hypomethylation have been demonstrated to play an important role in tumorigenesis.
Concentration:	1 mg/ml
Description:	Mouse monoclonal antibody to 5-methylcytosine (5-methylcytidine), clone 33D3.
Purification:	Protein A
Specificity:	Modified base 5-methylcytidine, a broad range of species.
Isotype:	IgG1
Formulation:	10 mM phosphate buffer, 150 mM NaCl, pH 7.4
Storage:	Aliquot and store at -20°C or lower. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance.
Application:	Dot Blot: 1:1000-1:2000; Immunohistochemistry: 1:100-1:500; Immunofluorescence: 1:100-1:500; ELISA: 1:1000-1:2000; MeDIP: 0.5-1 µg/reaction
Research Use:	Research use only.
Handling Recommendations:	For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Ordering Information:

Products	Size	Cat. No.
5-Methylcytosine (5-Methylcytidine)	10 µg	A-1014-010
Monoclonal Antibody [33D3]	50 µg	A-1014-050
	100 µg	A-1014-100

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