

DNMT1 Monoclonal Antibody [60B122.1]

(Catalog No. A-1001)

Background:	Dnmt1 has been demonstrated to preferably methylate hemimethylated DNA and is supposed to mediate both maintenance and <i>de novo</i> methylation. Increased activation or amount of Dnmt1 is believed to be involved in carcinogenesis, as well as other genetic and epigenetic diseases.
Concentration:	0.5 mg/ml
Description:	Mouse monoclonal antibody to Dnmt1, clone 60B1220.1 Epitope corresponding to amino acids 637-651 of human Dnmt1.
Purification:	Protein A Purified
Specificity:	Mouse and human.
Isotype:	IgG1
Formulation:	0.05% Sodium Azide, 0.2% Gelatin, PBS
Storage:	4°C, stable for 1 year from the date of shipment. For long-term storage, aliquot and store at -20°C.
Application:	Western Blot: 1:250 Immunofluorescence: 1:100 – 1:500 ELISA: 1:1000 – 1:2000 Immunoprecipitation: 2 µg/10 ⁶ cells
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.
DNMT1 Monoclonal Antibody [60B122.1]	50 µg	A-1001-050
	100 µg	A-1001-100
Available Related Products	Size	Cat. No.
DNMT3A Polyclonal Antibody	50 µg, 100 µg	A-1003
DNMT3B Polyclonal Antibody	50 µg, 100 µg	A-1004
DNMT3L Polyclonal Antibody	50 µg, 100 µg	A-1005

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