

MGMT Antibody [#C13084]

Catalog Number:	C13084
Concentration:	1mg/ml
Swiss-Prot No.:	P16455
Other Names:	Methylated-DNA--protein-cysteine methyltransferase, EC 2.1.1.63, 6-O-methylguanine-DNA methyltransferase, MGMT, O-6-methylguanine-DNA-alkyltransferase, MGMT
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human MGMT.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MGMT antibody detects endogenous levels of total MGMT protein.
Reactivity:	Human, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Tano K., Proc. Natl. Acad. Sci. U.S.A. 87:686-690(1990). Rydberg B., J. Biol. Chem. 265:9563-9569(1990). Koike G., J. Biol. Chem. 265:14754-14762(1990).