DNMT3B Antibody [#C11809]

Catalog Number: C11809
Concentration: 1mg/ml
Swiss-Prot No.: Q9UBC3

Other Names: DNA (cytosine-5)-methyltransferase 3B, EC 2.1.1.37, Dnmt3b, DNA methyltransferase HsallIB,

DNA MTase HsalliB, M.HsalliB, DNMT3B

Storage/Stability: Store at -20°C/1 year

Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02%

sodium azide and 50% glycerol.

Immunogen: The antiserum was produced against synthesized peptide derived from N-terminal of human

DNMT3B.

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-

specific immunogen.

Specificity: DNMT3B antibody detects endogenous levels of total DNMT3B protein.

Reactivity: Human

Applications: WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000

References: Deloukas P., Nature 414:865-871(2001).

Li H., J. Biol. Chem. 281:19489-19500(2006). Wijmenga C., Hum. Mutat. 16:509-517(2000).

Technical Support: support@aatbio.com; 408-733-1055