

NT5C3 Antibody [#C11674]

Catalog Number:	C11674
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H0P0
Other Names:	Cytosolic 5'-nucleotidase 3, EC 3.1.3.5, Cytosolic 5'-nucleotidase III, cN-III, Pyrimidine 5'-nucleotidase 1, P5'N-1, P5N-1, PN-I, Uridine 5'-monophosphate hydrolase 1, p36, NT5C3, P5N1, UMPH1, HSPC233
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 NT5C3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NT5C3 antibody detects endogenous levels of total NT5C3 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Amici A., Blood 96:1596-1598(2000). Marinaki A.M., Blood 97:3327-3332(2001). Kanno H., Br. J. Haematol. 126:265-271(2004).