

NDRG4 Antibody [#C18001]

Catalog Number:	C18001
Concentration:	1mg/ml
Swiss-Prot No.:	Q9ULP0
Other Names:	Brain development-related molecule 1; Protein NDRG4; SMAP-8; Vascular smooth muscle cell-associated protein 8
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 NDRG4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NDRG4 antibody detects endogenous levels of total NDRG4 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Zhou R.-H., Genomics 73:86-97(2001). Nishimoto S., Submitted (DEC-1998) to the EMBL/GenBank/DDBJ databases. Qu X., Mol. Cell. Biochem. 229:35-44(2002).