

NDRG3 Antibody [#C18000]

Catalog Number:	C18000
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UGV2
Other Names:	Protein NDRG3
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 NDRG3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NDRG3 antibody detects endogenous levels of total NDRG3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 92%); Rat (Identities = 92%, Positives = 92%)
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Zhou R.-H., Genomics 73:86-97(2001). Mao Y., Submitted (MAR-2000) to the EMBL/GenBank/DDBJ databases. Qu X., Mol. Cell. Biochem. 229:35-44(2002).