

NEDD8 Antibody [#C0359]

Catalog Number:	C0359
Concentration:	1mg/ml
Swiss-Prot No.:	Q15843
Other Names:	NED8, NEDD-8, Ubiquitin-like protein NEDD8
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 NEDD8.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NEDD8 antibody detects endogenous levels of total NEDD8 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Margaret A. Read, Mol. Cell. Biol., Apr 2000; 20: 2326 - 2333. Kenneth Wu, J. Biol. Chem., Aug 2003; 278: 28882 - 28891. Katsumi Kito, J. Biol. Chem., Jun 2001; 276: 20603 - 20609.