

SENP8 Antibody [#C0371]

Catalog Number:	C0371
Concentration:	1mg/ml
Swiss-Prot No.:	Q96LD8
Other Names:	DEN1, NEDP, DENEDDYLASE 1, NEDD8-SPECIFIC PROTEASE 1, SENTRIN-SPECIFIC PROTEASE FAMILY, MEMBER 8, SUMO/sentrin specific protease family member 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 SENP8.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SENP8 antibody detects endogenous levels of total SENP8 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	Tudeviin Gan-Erdene, J. Biol. Chem., Aug 2003; 278: 28892 - 28900. Debaditya Mukhopadhyay, J. Cell Biol., Sep 2006; 174: 939 - 949. Thomas Colby, Plant Physiology, Sep 2006; 142: 318 - 332. Joris Hemelaar, Mol. Cell. Biol., Jan 2004; 24: 84 - 95.