

AOS1 Antibody [#C0351]

Catalog Number:	C0351
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UBE0
Other Names:	AOS1, HSPC140, SAE1, SUA1, SUMO-1 activating enzyme E1 N subunit, ULE1A, UBLE1A
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 AOS1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AOS1 antibody detects endogenous levels of total AOS1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Yoshiaki Azuma, FASEB J, Jun 2001; 10.1096. Walter Makous, Invest. Ophthalmol. Vis. Sci., Sep 2006; 47: 4160 - 4167. Hsin-Pei Shih, J. Cell Sci., Mar 2002; 115: 1259. Thomas Buschmann, J. Biol. Chem., Oct 2001; 276: 40389 - 40395.