

COPS5 Antibody [#C11442]

Catalog Number:	C11442
Concentration:	1mg/ml
Swiss-Prot No.:	Q92905
Other Names:	EC 3.4.-.-, Signalosome subunit 5, SGN5, Jun activation domain-binding protein 1
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 COPS5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	COPS5 antibody detects endogenous levels of total COPS5 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Claret F.-X., Nature 383:453-457(1996). Asano K., J. Biol. Chem. 272:27042-27052(1997). The MGC Project Team, Genome Res. 14:2121-2127(2004).