

GPSN2 Antibody [#C18998]

Catalog Number:	C18998
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NZ01
Other Names:	Synaptic glycoprotein SC2; EC 1.3.1.-; Trans-2,3-enoyl-CoA reductase; TER; GPSN2; SC2
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 GPSN2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPSN2 antibody detects endogenous levels of total GPSN2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Mao M., Proc. Natl. Acad. Sci. U.S.A. 95:8175-8180(1998). Seo H.C., Submitted (JAN-2000) to the EMBL/GenBank/DDBJ databases. Kalnine N., Submitted (MAY-2003) to the EMBL/GenBank/DDBJ databases.