

MYOM1 Antibody [#C16759]

Catalog Number:	C16759
Concentration:	1mg/ml
Swiss-Prot No.:	P52179
Other Names:	Myomesin-1, Myomesin family member 1; 190 kDa titin-associated protein; 190 kDa connectin-associated protein
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 MYOM1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MYOM1 antibody detects endogenous levels of total MYOM1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Hayess K., Submitted (JAN-2004) to the EMBL/GenBank/DDBJ databases. Ota T., Nat. Genet. 36:40-45(2004). Nusbaum C., Nature 437:551-555(2005).