

MYOM2 Antibody [#C16760]

Catalog Number:	C16760
Concentration:	1mg/ml
Swiss-Prot No.:	P54296
Other Names:	165 kDa connectin-associated protein; 165 kDa titin-associated protein; M-protein; MYOM2; myomesin-2
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 MYOM2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MYOM2 antibody detects endogenous levels of total MYOM2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Vinkemeier U., J. Cell Sci. 106:319-330(1993).