

## MYSM1 Antibody [#C17997]

<b>Catalog Number:</b>	C17997
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q5VVJ2
<b>Other Names:</b>	Myb-like, SWIRM and MPN domain-containing protein 1
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human MYSM1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	MYSM1 antibody detects endogenous levels of total MYSM1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100   ELISA: 1:40000
<b>References:</b>	Nagase T., DNA Res. 8:179-187(2001). Gregory S.G., Nature 441:315-321(2006).Matsuoka S.,Science 316:1160-1166(2007).