

JMY Antibody [#C10003]

Catalog Number:	C10003
Concentration:	1mg/ml
Swiss-Prot No.:	A1L4P5
Other Names:	JMY protein, JMY
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 JMY.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	JMY antibody detects endogenous levels of total JMY protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Strausberg R.L., Proc. Natl. Acad. Sci. U.S.A. 99:16899-16903(2002). NIH MGC Project, Submitted (JAN-2007) to the EMBL/GenBank/DDBJ databases.