

TMBIM1 Antibody [#C19281]

Catalog Number:	C19281
Concentration:	1mg/ml
Swiss-Prot No.:	Q969X1
Other Names:	PP1201; PSEC0158; RECS1; TMBI1; transmembrane BAX inhibitor motif containing 1; transmembrane BAX inhibitor motif-containing 1
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 TMBIM1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TMBIM1 antibody detects endogenous levels of total TMBIM1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Ota T., Submitted (MAR-2002) to the EMBL/GenBank/DDBJ databases. Wan D., Proc. Natl. Acad. Sci. U.S.A. 101:15724-15729(2004). Ota T., Nat. Genet. 36:40-45(2004).