

MYST1 Antibody [#C11704]

Catalog Number:	C11704
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H7Z6
Other Names:	EC 2.3.1.48; MYST protein 1; MOZ, YBF2/SAS3, SAS2 and TIP60 protein 1; hMOF
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 MYST1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MYST1 antibody detects endogenous levels of total MYST1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	Wan D., Proc. Natl. Acad. Sci. U.S.A. 101:15724-15729(2004). Neal K.C., Biochim. Biophys. Acta 1490:170-174(2000). Pardo P.S., J. Biol. Chem. 277:50860-50866(2002).