

## MYST2 Antibody [#C10221]

<b>Catalog Number:</b>	C10221
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O95251
<b>Other Names:</b>	EC 2.3.1.48, HBO1, HBOA, MYST histone acetyltransferase 2
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), 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 MYST2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	MYST2 antibody detects endogenous levels of total MYST2 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Iizuka M., J. Biol. Chem. 274:23027-23034(1999). Sharma M., J. Biol. Chem. 275:35200-35208(2000). Olsen J.V., Cell 127:635-648(2006).