

B-Myb (Ab-577/581) Antibody [#B0823]

Catalog Number:	B0823
Concentration:	1mg/ml
Swiss-Prot No.:	P10244
Other Names:	BMYP, MYBB, MYBL2, Myb-related protein B
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 non-phosphopeptide derived from human B-Myb around the phosphorylation site of serine 577/581 (R-R-S ^P -P-I).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	B-Myb (Ab-577/581) antibody detects endogenous levels of total B-Myb protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Nomura N., Nucleic Acids Res. 16:11075-11089(1988). Deloukas P., Nature 414:865-871(2001). Olsen J.V., Cell 127:635-648(2006). Litovchick L., Mol. Cell 26:539-551(2007).