

Myc Antibody [#C10262]

Catalog Number:	C10262
Concentration:	1mg/ml
Swiss-Prot No.:	P01106
Other Names:	Myc proto-oncogene protein, c-myc
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 C-terminal of human Myc.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Myc antibody detects endogenous levels of total Myc protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Batley J., Cell 34:779-787(1983). Bernard O., EMBO J. 2:2375-2383(1983). Colby W.W., Nature 301:722-725(1983).