

Myc (Ab-58) Antibody [#B7159]

- Catalog Number:** B7159-1, B7159-2, B7159
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- 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.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human Myc around the phosphorylation site of threonine 58.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** Myc (Ab-58) antibody detects endogenous levels of total Myc protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:20000
- Swiss-Prot No. :** P01106
- References:** Jin Z, et al. (2004) J Biol Chem. 279(38): 40209-40219.
Welcker M, et al. (2004) Proc Natl Acad Sci U S A. 101(24): 9085-9090.
Baudino T A, et al. (2001) Mol Cell Biol. 21: 691-702.
Blackwood E M, et al. (1991) Science. 251:1211-1217.
Henriksson M, et al. (1996) Adv Cancer Res. 68: 109-182.