

IL3R (Ab-593) Antibody [#B0444]

- Catalog Number:** B0444-1, B0444-2, B0444
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P26951
- All Names:** Interleukin-3 receptor alpha chain, IL-3R-alpha, CD123 antigen, IL3RA
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), 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 IL3R around the phosphorylation site of tyrosine 593 (G-P-Y-L-G).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** IL3R (Ab-593) antibody detects endogenous levels of total IL3R protein.
- Reactivity:** Human, Mouse
- Applications:** WB: 1:500~1:1000 ELISA: 1:5000
- References:** Michelle M. Le Beau, Blood, Aug 2003; 102: 1072 - 1074.
Elena Cattaneo, J. Biol. Chem., Sep 1996; 271: 23374.
Patricia McLoughlin, J. Biol. Chem., Sep 2002; 277: 37045 - 37053.
Louise Edvardsson, Blood, Nov 2004; 104: 4120.