

Actin- α -1 Antibody [#C0121]

- Catalog Number:** C0121
- Amount:** 50 μ g/50 μ l, 100 μ g/100 μ l, 200 μ g/200 μ l
- Swiss-Prot No. :** P68133
- All Names:** ACTA, ACTA1, ACTS, Actin, alpha skeletal muscle, Alpha-actin 1
- 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 peptide derived from human Actin- α -1.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Actin- α -1 antibody detects endogenous levels of total Actin- α -1 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
- References:** Atsushi Masamune, J. Pharmacol. Exp. Ther., Jan 2003; 304
Alicia Algeciras-Schimmich, PNAS, Sep 2003; 100: 11445.
Richard D. Peavy, Mol. Pharmacol., Jun 2005; 67: 2102 - 2114.
Xiaobing Deng, J. Biol. Chem., Feb 2005; 280: 4894 - 4905.