

Actin-pan ($\alpha/\beta/\gamma$) (Ab-53/55) Antibody [#B0762]

Catalog Number:	B0762
Concentration:	1mg/ml
Swiss-Prot No.:	P60709/P63261
Other Names:	ACTA, ACTA1, ACTS, Actin, alpha skeletal muscle, Alpha-actin 1; ACTB, Actin, cytoplasmic 1, Beta-actin; ACTG, ACTG1, actin, cytoplasmic 2, gamma-actin
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 Actin-pan ($\alpha/\beta/\gamma$) around the phosphorylation site of tyrosine 53/55 (D-S-Y ^P -V-G).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Actin-pan ($\alpha/\beta/\gamma$) (Ab-53/55) antibody detects endogenous levels of total Actin-pan ($\alpha/\beta/\gamma$) protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	A. J. MacQueen, Mol. Biol. Cell, Jul 2005; 16: 3247 - 3259. Isaac Rabinovitz, J. Cell Biol., Dec 1997; 139: 1873. Deanne H. Hryciw, Am J Physiol Renal Physiol, Jun 2005; 288: F1227 - F1235. Derek T. Warren, J. Cell Sci., Apr 2002; 115: 1703.