

YAP (Ab-127) Antibody [#B0757]

Catalog Number:	B0757
Concentration:	1mg/ml
Swiss-Prot No.:	P46937
Other Names:	65 kDa Yes-associated protein, YAP1, YAP65
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 YAP around the phosphorylation site of serine 127 (A-H-S ^P -S-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	YAP (Ab-127) antibody detects endogenous levels of total YAP protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Rami I. Aqeilan, Cancer Res., Aug 2005; 65: 6764 - 6772. Michael Howell, J. Biol. Chem., Jun 2004; 279: 26300 - 26306. Akihiko Komuro, J. Biol. Chem., Aug 2003; 278: 33334 - 33341. Masataka Hoshino, J. Cell Biol., Feb 2006; 172: 589 - 604.