

mTOR (Ab-2446) Antibody [#B1156]

Catalog Number:	B1156
Concentration:	1mg/ml
Swiss-Prot No.:	P42345
Other Names:	FRAP, FRAP1, FRAP2, RAFT1, Rapamycin target protein, kinase mTOR
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 mTOR around the phosphorylation site of threonine 2446 (T-R-T ^P -D-S).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	mTOR (Ab-2446) antibody detects endogenous levels of total mTOR protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Brown E.J., Nature 369:756-758(1994). Onyango P., Genomics 50:187-198(1998). Gregory S.G., Nature 441:315-321(2006).