

Tuberin (Ab-981) Antibody [#B8443]

Catalog Number:	B8443
Concentration:	1mg/ml
Swiss-Prot No.:	P49815
Other Names:	TSC2; tuberous sclerosis 2 gene; Tuberous sclerosis 2 homolog protein; Tuberous sclerosis 2 protein
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 Tuberin around the phosphorylation site of serine 981 (S-I-S ^P -V-S).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Tuberin (Ab-981) antibody detects endogenous levels of total Tuberin protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Nellist M., Cell 75:1305-1315(1993). Xu L., Genomics 27:475-480(1995). Martin J., Nature 432:988-994(2004).