

Tuberin/TSC2 Antibody [#B0591]

Catalog Number:	B0591
Concentration:	1mg/ml
Swiss-Prot No.:	P49815
Other Names:	TSC2, 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 peptide derived from Internal of human Tuberin/TSC2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Tuberin/TSC2 antibody detects endogenous levels of total Tuberin/TSC2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
References:	Ralf Wienecke, J. Biol. Chem., Jul 1995; 270: 16409. Lauri D. Aicher, J. Biol. Chem., Jun 2001; 276: 21017 - 21021. Geraldine A. Finlay, J. Biol. Chem., May 2004; 279: 23114 - 23122. Vanishree Murthy, Am J Physiol Renal Physiol, May 2000; 278: 737.