

ITCH (Ab-420) Antibody [#B8254]

Catalog Number:	B8254
Concentration:	1mg/ml
Swiss-Prot No.:	Q96J02
Other Names:	AIF4; AIP4; atrophin-1-interacting protein 4; EC 6.3.2.-; Itchy homolog E3 ubiquitin protein ligase; NAPP1; NFE2-associated polypeptide 1; ubiquitin protein ligase ITCH
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 ITCH around the phosphorylation site of tyrosine 420 (F-I-Y ^P -G-N).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ITCH (Ab-420) antibody detects endogenous levels of total ITCH protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Chen X., Genomics 73:238-241(2001). Deloukas P., Nature 414:865-871(2001). Wood J.D., Mol. Cell. Neurosci. 11:149-160(1998).