

TLK1 Antibody [#B0749]

Catalog Number:	B0749
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y4F6
Other Names:	EC 2.7.11.1, PKU-beta, TOUSLED-like kinase 1 homolog, Tousled-like kinase 1, kinase TLK1
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 TLK1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TLK1 antibody detects endogenous levels of total TLK1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Yamakawa A., Gene 202:193-201(1997). Sillje H.H.W., EMBO J. 18:5691-5702(1999). Nakajima D., DNA Res. 9:99-106(2002)