

MRCKB Antibody [#C11973]

Catalog Number:	C11973
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y5S2
Other Names:	CDC42-binding protein kinase beta, CDC42BPB, EC 2.7.11.1, MRCK-beta
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 MRCKB.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRCKB antibody detects endogenous levels of total MRCKB protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Nakajima D., DNA Res. 9:99-106(2002). Olsen J.V., Cell 127:635-648(2006). Hirosawa M., DNA Res. 6:329-336(1999).