

KV2.1/KCNB1 Antibody [#B1086]

Catalog Number:	B1086
Concentration:	1mg/ml
Swiss-Prot No.:	Q14721
Other Names:	DRK1, KCB1, KCNB1, Potassium channel Kv2.1, Potassium voltage-gated channel subfamily B member 1
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 KV2.1/KCNB1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KV2.1/KCNB1 antibody detects endogenous levels of total KV2.1/KCNB1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 90%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	G Shi, J. Biol. Chem., Sep 1994; 269: 23204 - 23211. Jobst-Hendrik Schultz, Circ. Res., Mar 2001; 88: 483 - 490. RH Scannevin, J. Cell Biol., Dec 1996; 135: 1619.