

KCNV2 Antibody [#C17829]

Catalog Number:	C17829
Concentration:	1mg/ml
Swiss-Prot No.:	Q8TDN2
Other Names:	KCNV2; potassium voltage-gated channel subfamily V member 2; voltage-gated potassium channel subunit Kv8.2
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 KCNV2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KCNV2 antibody detects endogenous levels of total KCNV2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 87%, Positives = 93%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Ottshytsch N., Proc. Natl. Acad. Sci. U.S.A. 99:7986-7991(2002). Humphray S.J., Nature 429:369-374(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004).