

## Kv1.3/KCNA3 Antibody [#B1085]

<b>Catalog Number:</b>	B1085
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P22001
<b>Other Names:</b>	CIK3, HGK5, HLK3, HPCN3, HuKIII, KCNA3, Potassium channel Kv1.3, Potassium voltage-gated channel subfamily A member 3
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from Internal of human Kv1.3/KCNA3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Kv1.3/KCNA3 antibody detects endogenous levels of total Kv1.3/KCNA3 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:100~1:500 ELISA: 1:20000
<b>References:</b>	Rubén Vicente, J. Biol. Chem., Dec 2006; 281: 37675 - 37685. Christine Beeton, PNAS, Nov 2006; 103: 17414 - 17419. Christopher B. Fordyce, J. Neurosci., Aug 2005; 25: 7139 - 7149.