

GRK2 Antibody [#B0486]

Catalog Number:	B0486
Concentration:	1mg/ml
Swiss-Prot No.:	P25098
Other Names:	ADRBK1, ARBK1, BARK, BARK1, Beta-ARK-1, Beta-adrenergic receptor kinase 1, EC 2.7.11.15, G protein receptor kinase 2, G- protein coupled receptor kinase 2, kinase GRK2
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 GRK2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GRK2 antibody detects endogenous levels of total GRK2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
References:	Aarti N. Desai, J. Pharmacol. Exp. Ther., Jun 2006; 317: 1027 - 1035. John A. Gray, Mol. Pharmacol., Nov 2001; 60: 1020. Anne Vroon, J. Leukoc. Biol., May 2004; 75: 901 - 909. R. J. Hughes, Am J Physiol Cell Physiol, Mar 1996; 270: C885 - C891.