

DCLK3 Antibody [#C11659]

Catalog Number :	C11659
Concentration:	1mg/ml
Swiss-Prot No.:	Q9C098
Other Names:	Serine/threonine-protein kinase DCLK3, EC 2.7.11.1, Doublecortin-like and CAM kinase-like 3, Doublecortin-like kinase 3 , DCLK3, DCAMKL3, DCDC3C, KIAA1765
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 DCLK3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	DCLK3 antibody detects endogenous levels of total DCLK3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
References:	Nagase T., DNA Res. 7:347-355(2000). Ota T., Nat. Genet. 36:40-45(2004).