

Hexokinase-3 Antibody [#C10708]

Catalog Number:	C10708
Concentration:	1mg/ml
Swiss-Prot No.:	P52790
Other Names:	Hexokinase-3, EC 2.7.1.1, Hexokinase type III, HK III, HK3
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 Hexokinase-3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Hexokinase-3 antibody detects endogenous levels of total Hexokinase-3 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Furuta H., Genomics 36:206-209(1996). Palma F., Mol. Cell. Biochem. 155:23-29(1996). Sjoebloom T., Science 314:268-274(2006).