

NBPF5 Antibody [#C16930]

Catalog Number:	C16930
Concentration:	1mg/ml
Swiss-Prot No.:	Q86XG9
Other Names:	Neuroblastoma breakpoint family member 5
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 C-terminal of human NBPF5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NBPF5 antibody detects endogenous levels of total NBPF5 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	The MGC Project Team; Genome Res. 14-2121-2127(2004). Vandepoele K., Mol. Biol. Evol. 22:2265-2274(2005).