

PHF3 Antibody [#C17658]

Catalog Number:	C17658
Concentration:	1mg/ml
Swiss-Prot No.:	Q92576
Other Names:	PHD finger protein 3
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 PHF3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PHF3 antibody detects endogenous levels of total PHF3 protein.
Reactivity:	Human (Identities = 92%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Fischer U., Cytogenet. Cell Genet. 94:131-136(2001). Nagase T., DNA Res. 3:321-329(1996). Nakajima D., DNA Res. 9:99-106(2002).