

PHLA1 Antibody [#C13101]

Catalog Number:	C13101
Concentration:	1mg/ml
Swiss-Prot No.:	Q8WV24
Other Names:	Pleckstrin homology-like domain family A member 1; T-cell death-associated gene 51 protein; Apoptosis-associated nuclear protein; Proline- and histidine-rich protein; Proline- and glutamine-rich protein; PQ-rich protein; PHLDA1; PHRIP, TDAG51
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 PHLA1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PHLA1 antibody detects endogenous levels of total PHLA1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Kuske M.D., Cytogenet. Cell Genet. 89:1-1(2000). Neef R., Cancer Res. 62:5920-5929(2002). Hinz T., Cell. Signal. 13:345-352(2001).