

PARP4 Antibody [#C13099]

Catalog Number:	C13099
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UKK3
Other Names:	193-kDa vault protein, ADP- ribose, ADP-ribosyltransferase (NAD ⁺ , ADPRTL1, EC 2.4.2.30, H5 proline-rich, I-alpha-I-related, PARP- related/lalphal-related H5/proline-rich, PARP-4, PARP-related, PARPL, PH5P, VAULT3, VPARP, p193, poly (ADP-ribose) polymerase family, member 4, poly (ADP-ribose) polymerase)-like 1, poly [ADP-ribose] polymerase 4, poly(ADP-ribose) synthetase, poly(ADP-ribosyl)transferase-like 1, polymerase, vault poly, vault protein, 193-kDa
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 PARP4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PARP4 antibody detects endogenous levels of total PARP4 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	Kickhoefer V.A., J. Cell Biol. 146:917-928(1999). Dunham A., Nature 428:522-528(2004). Still I.H., Genomics 62:533-536(1999).