

TP53INP2 Antibody [#C11320]

Catalog Number:	C11320
Concentration:	1mg/ml
Swiss-Prot No.:	Q8CFU8
Other Names:	p53-inducible protein U; PIG-U; PINH; T53I2; tumor protein p53 inducible nuclear protein 2; tumor protein p53-inducible nuclear protein 2
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 TP53INP2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TP53INP2 antibody detects endogenous levels of total TP53INP2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Baumgartner B.G., Submitted (OCT-2000) to the EMBL/GenBank/DDBJ databases. Deloukas P., Nature 414:865-871(2001). The MGC Project Team, Genome Res. 14:2121-2127(2004).