

TP53INP1 Antibody [#C11471]

Catalog Number:	C11471
Concentration:	1mg/ml
Swiss-Prot No.:	Q96A56
Other Names:	Tumor protein p53-inducible nuclear protein 1, p53-dependent damage-inducible nuclear protein 1, p53DINP1, Stress-induced protein, TP53INP1, P53DINP1, SIP
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 TP53INP1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TP53INP1 antibody detects endogenous levels of total TP53INP1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Okamura S., Mol. Cell 8:85-94(2001). Tomasini R., Eur. J. Cell Biol. 81:294-301(2002). Tomasini R., J. Biol. Chem. 278:37722-37729(2003).