

WWOX Antibody [#C13129]

Catalog Number:	C13129
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NZC7
Other Names:	WW domain-containing oxidoreductase, EC 1.1.1.-, Fragile site FRA16D oxidoreductase, WWOX, FOR, WOX1
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 WWOX.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	WWOX antibody detects endogenous levels of total WWOX protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Bednarek A.K., Cancer Res. 60:2140-2145(2000). Ried K., Hum. Mol. Genet. 9:1651-1663(2000). Bednarek A.K., Cancer Res. 61:8068-8073(2001).