

OXR1 Antibody [#C17579]

Catalog Number:	C17579
Concentration:	1mg/ml
Swiss-Prot No.:	Q8N573
Other Names:	Oxidation resistance protein 1
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 OXR1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	OXR1 antibody detects endogenous levels of total OXR1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Ohira M., Cancer Lett. 197:63-68(2003). The German cDNA consortium; Submitted (JUN-2005) to the EMBL/GenBank/DDBJ databases.