

MOX2R Antibody [#C20886]

Catalog Number:	C20886
Concentration:	1mg/ml
Swiss-Prot No.:	Q8TD46
Other Names:	CD200 cell surface glycoprotein receptor; CD200R1; Cell surface glycoprotein OX2 receptor precursor; CRTR2; MOX2R; OX2R
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 MOX2R.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MOX2R antibody detects endogenous levels of total MOX2R protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Gorczynski R.M., Am. J. Reprod. Immunol. 52:147-163(2004). Wright G.J., Submitted (JUN-2000) to the EMBL/GenBank/DDBJ databases. Suarez A., Submitted (APR-2002) to the EMBL/GenBank/DDBJ databases.