

Aquaporin 2 Antibody [#B0768]

Catalog Number:	B0768
Concentration:	1mg/ml
Swiss-Prot No.:	P41181
Other Names:	ADH water channel, AQP-CD, Aquaporin-CD, Collecting duct water channel protein, WCH-CD, Water channel protein for renal collecting duct, aquaporin 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 Aquaporin 2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Aquaporin 2 antibody detects endogenous levels of total Aquaporin 2 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: 1:10000
References:	Sasaki S., J. Clin. Invest. 93:1250-1256(1994). van Balkom B.W.M., J. Biol. Chem. 277:41473-41479(2002). Oksche A., Hum. Genet. 98:587-589(1996).