

Aquaporin 2(Phospho-Ser256)Antibody

Catalog Number:	8A0768
Concentration:	1mg/ml
Origin of Antibody:	Rabbit
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
All Sites:	Human: Ser256; Mouse: Ser256; Rat: Ser256
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 phosphopeptide derived from human Aquaporin 2 around the phosphorylation site of serine 256 (R-Q-S ^P -V-E).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity:	Aquaporin 2 (Phospho-Ser256) antibody detects endogenous levels of Aquaporin 2 only when phosphorylated at serine 256.
Reactivity:	Human, Mouse, Rat.
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Takeaki Inoue, Am J Physiol Renal Physiol, Nov 1998; 275: 752. PM Deen, J. Am. Soc. Nephrol., Oct 1997; 8: 1493. Baoxue Yang, J. Biol. Chem., Jan 2001; 276: 2775 - 2779.