

## AQP1 Antibody [#C14546]

<b>Catalog Number:</b>	C14546
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P29972
<b>Other Names:</b>	AQP-1; AQP-CHIP; aquaporin 1; aquaporin-1; aquaporin-CHIP; channel-like integral membrane protein, 28-kDa; CHIP28; Colton blood group; urine water channel; water channel protein for red blood cells and kidney proximal tubule
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human AQP1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	AQP1 antibody detects endogenous levels of total AQP1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 92%); Rat (Identities = 92%, Positives = 92%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Preston G.M., Proc. Natl. Acad. Sci. U.S.A. 88:11110-11114(1991). Moon C., J. Biol. Chem. 268:15772-15778(1993). Ruiz A.C., Biochim. Biophys. Acta 1282:174-178(1996).