

## Claudin 1 (Ab-210) Antibody [#B8316]

<b>Catalog Number:</b>	B8316
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O95832
<b>Other Names:</b>	CLD1; CLDN1; SEMP1; Senescence-associated epithelial membrane protein; UNQ481/PRO944
<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 non-phosphopeptide derived from human Claudin 1 around the phosphorylation site of tyrosine 210 (G-K-D-Y <sup>P</sup> -V).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Claudin 1 (Ab-210) antibody detects endogenous levels of total Claudin 1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Swisshelm K.L., Gene 226:285-295(1999). Halford S., Cytogenet. Cell Genet. 88:217-217(2000). Kraemer F., Hum. Genet. 107:249-256(2000).