

Claudin 3 (Ab-219) Antibody [#B8317]

Catalog Number:	B8317
Concentration:	1mg/ml
Swiss-Prot No.:	O15551
Other Names:	CLD3; CLDN3; Clostridium perfringens enterotoxin receptor 2; CPE-R 2; CPE-receptor 2; HRVP1; Ventral prostate 1 protein homolog
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 non-phosphopeptide derived from human Claudin 3 around the phosphorylation site of tyrosine 219 (R-K-D-Y ^P -V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Claudin 3 (Ab-219) antibody detects endogenous levels of total Claudin 3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Peacock R.E., Genomics 46:443-449(1997). Katahira J., J. Biol. Chem. 272:26652-26658(1997). The MGC Project Team; Genome Res. 14:2121-2127(2004).