

Claudin 1 Antibody [#C0142]

Catalog Number: C0142-1, C0142-2, C0142

Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl

Swiss-Prot No. : O95832

All Names: CLD1, CLDN1, SEMP1, Senescence-associated epithelial membrane protein

Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Storage/Stability: Store at -20°C/1 year

Immunogen: The antiserum was produced against synthesized peptide derived from human Claudin 1.

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.

Specificity/Sensitivity Claudin 1 antibody detects endogenous levels of total Claudin 1 protein.

Reactivity: Human, Mouse, Rat

Applications: WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000

References: Julie Dufresne and Daniel G. Cyr, Biol Reprod, May 2007; 76: 825 - 832.

Evemie Dubé, Biol Reprod, Apr 2007; 10.1095/biolreprod.106.059246.

Michele A. Battle, PNAS, May 2006; 103: 8419 - 8424.

Edmund C. Chang, Endocrinology, Oct 2006; 147: 4831 - 4842.