

E-cadherin-pan Antibody [#C0148]

- Catalog Number:** C0148-1, C0148-2, C0148
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P12830
- All Names:** CADH1, CAM 120/80, CDH1, cadherin-1, epithelial-cadherin, uvomorulin
- 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 E-cadherin.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** E-cadherin-pan antibody detects endogenous levels of total E-cadherin protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
- References:** Michaël Perrais, Mol. Biol. Cell, Mar 2007; 10.1091/mbc.E06-04-0348.
Fiorella Balzac, J. Cell Sci., Oct 2005; 118: 4765 - 4783.
Gianpaolo Suriano, Hum. Mol. Genet., Nov 2003; 12: 3007 - 3016.
Seiyu Hosono, J. Biol. Chem., Apr 2000; 275: 10943.