

Catenin- α Antibody [#C0137]

- Catalog Number:** C0137
- Amount:** 50 μ g/50 μ l, 100 μ g/100 μ l, 200 μ g/200 μ l
- Swiss-Prot No. :** P35221
- All Names:** 102 kDa cadherin-associated protein, Alpha E-catenin, CATNA1, CTN1, Cadherin-associated protein, catenin alpha-1
- 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 Catenin- α .
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Catenin- α antibody detects endogenous levels of total Catenin- α protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
- References:** Yutaka Shimoyama, Cancer Res., Oct 1992; 52: 5770 - 5774.
Griet Vanpoucke, Nucleic Acids Res., Aug 2004; 32: 4155 - 4165.
Jolanda van Hengel, J. Cell Biol., Jun 1997; 137: 1103.
Masahiko Itoh, J. Cell Biol., Jul 1997; 138: 181.