

Keratin 10 Antibody [#C0249]

- Catalog Number:** C0249-1, C0249-2, C0249
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P13645
- All Names:** Keratin type I cytoskeletal 10, Cytokeratin-10, CK-10, K10
- 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 Keratin 10.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Keratin 10 antibody detects endogenous levels of total Keratin 10 protein.
- Reactivity:** Human, Rat
- Applications:** WB: 1:300~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
- References:** XM Zhou, J. Biol. Chem., Oct 1988; 263: 15584 - 15589.
Felix B. Müller, Hum. Mol. Genet., Apr 2006; 15: 1133 - 1141.
J Reichelt, J. Cell Sci., Sep 1997; 110: 2175.
BP Korge, PNAS, Feb 1992; 89: 910.