

Keratin 7 Antibody [#C0247]

- Catalog Number:** C0247-1, C0247-2, C0247
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P08729
- All Names:** Keratin type II cytoskeletal 7, Cytokeratin-7, CK-7, K7, Sarcolectin
- 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 7.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Keratin 7 antibody detects endogenous levels of total Keratin 7 protein.
- Reactivity:** Human
- Applications:** WB: 1:300~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
- References:** M Rosenberg, Mol. Cell. Biol., Feb 1988; 8: 722 - 736.
SM Troyanovsky, J. Cell Sci., Jul 1989; 93: 419.
Kevin M. McGowan, Genes & Dev., Jun 2002; 16: 1412.
Nicholas Bianchi, Mol. Cell. Biol., Aug 2005; 25: 7249 - 7259.