

Keratin 19 Antibody [#C0244]

Catalog Number: C0244-1, C0244-2, C0244

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

Swiss-Prot No. : P08727

All Names: Keratin type I cytoskeletal 19, Cytokeratin-19, CK-19, K19

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 19.

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

Specificity/Sensitivity Keratin 19 antibody detects endogenous levels of total Keratin 19 protein.

Reactivity: Human, Mouse, Rat

Applications: WB: 1:300~1:1000 IHC: 1:50~1:100 ELISA: 1:10000

References: Lianna Wong, Chem Senses, Jan 1994; 19: 251 - 264.

L Hu and LJ Gudas J. Biol. Chem., Jan 1994; 269: 183 - 191.

Yong-Sik Jung, Jpn. J. Clin. Oncol., Apr 2003; 33: 167 - 172.

Xiangjun Zhou, J. Biol. Chem., Apr 1999; 274: 12861.