

## p63 Antibody [#C0292]

**Catalog Number:** C0292-1, C0292-2, C0292

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

**Swiss-Prot No. :** Q9H3D4

**All Names:** EEC3, KET, LMS, SHFM4, TA p63 alpha, TP63, p51, p73H, p73L, tumor protein p63

**Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg<sup>2+</sup> and Ca<sup>2+</sup>), 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 p63 antibody.

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

**Specificity/Sensitivity** p63 antibody detects endogenous levels of total p63 protein.

**Reactivity:** Human, Mouse, Rat

**Applications:** WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000

**References:** Pranay C. Patel, Arch Otolaryngol Head Neck Surg, May 1998; 124: 559.

Atul A. Gawande, Arch Surg, Apr 2006; 141: 381 - 384.

Timothy M. Murray, Endocr. Rev., Feb 2005; 26: 78 - 113.

W. M. Philbrick, Physiol Rev, Jan 1996; 76: 127.