

## p15 INK Antibody [#C0287]

**Catalog Number:** C0287-1, C0287-2, C0287

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

**Swiss-Prot No. :** P42772

**All Names:** CDKN2B, CDN2B, Cyclin-dependent kinase 4 inhibitor B, MTS2, Multiple tumor suppressor 2, p14-INK4b

**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 p15 INK antibody.

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

**Specificity/Sensitivity** p15 INK antibody detects endogenous levels of total p15 INK protein.

**Reactivity:** Human, Mouse, Rat

**Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000

**References:** Nynke Y. Rots, Blood, Apr 1999; 93: 2721 - 2729.

Stefan Faderl, Clin. Cancer Res., Jul 1999; 5: 1855.

George Wu, Mol. Biol. Cell, Mar 2007; 18: 1018 - 1029.

A. Koenig, Vet. Pathol., Jul 2002; 39: 458.