

p18 INK Antibody [#C0288]

Catalog Number: C0288-1, C0288-2, C0288

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

Swiss-Prot No. : P42773

All Names: OTTHUMP00000046546, p18(INK4C), p18-INK4C, CDK6 inhibitor p18, cyclin-dependent inhibitor, cyclin-dependent kinase 4 inhibitor C, cyclin-dependent kinase 6 inhibitor p18, cyclin-dependent kinase inhibitor 2C (p18, inhibits CDK4)

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 p18 INK antibody.

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

Specificity/Sensitivity p18 INK antibody detects endogenous levels of total p18 INK protein.

Reactivity: Human, Mouse, Rat

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

References: Youzhong Yuan, Cancer Res., Jan 2006; 66: 343 - 351.

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

Pei-Fang Hung, Am J Physiol Cell Physiol, May 2005; 288: C1094 - C1108.

Mana Mirjany, J. Pharmacol. Exp. Ther., May 2002; 301: 494.