

## Rb (Ab-795) Antibody [#B7209]

- Catalog Number:** B7209-1, B7209-2, B7209
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- 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 non-phosphopeptide derived from human Rb around the phosphorylation site of serine 795 (P-S-S<sup>P</sup>-P-L).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** Rb (Ab-795) antibody detects endogenous levels of total Rb protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000    ELISA: 1:20000
- Swiss-Prot No. :** P06400
- References:** Roesch A, et al. (2005) *Mod Pathol.* 18(4): 565-572.  
Chadee DN, et al. (2004) *Nat Cell Biol.* 6(8): 770-776.  
Knudsen ES, et al. (1997) *Mol Cell Biol.* 17(10): 5771-5783.  
Knudsen ES, et al. (1996) *J Biol Chem.* 271(14): 8313-8320.