

RIT1 Antibody [#C0317]

- Catalog Number:** C0317-1, C0317-2, C0317
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** Q92963
- All Names:** GTP-binding protein Rit1, Ras-like protein expressed in many tissues, Ras-like without CAAX protein 1, RIBB, RIT, ROC1
- 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 RIT1.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** RIT1 antibody detects endogenous levels of total RIT1 protein.
- Reactivity:** Human, Mouse
- Applications:** WB: 1:500~1:1000 ELISA: 1:10000
- References:** Manabu Furukawa, Mol. Cell. Biol., Nov 2000; 20: 8185 - 8197.
Angus Chen, J. Biol. Chem., May 2000; 275: 15432.
Kenneth Wu, Mol. Cell. Biol., Feb 2000; 20: 1382 - 1393.
Kenneth Wu, J. Biol. Chem., Jan 2002; 277: 516 - 527.