

RAN Antibody [#C0312]

- Catalog Number:** C0312-1, C0312-2, C0312
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Swiss-Prot No. : P62826
All Names: ARA24, Androgen receptor- associated protein 24, LPS, RASL2-8, Ran GTPase, TC4
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 RAN.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity RAN antibody detects endogenous levels of total RAN protein.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 ELISA: 1:10000
References: Karen M. Lounsbury, J. Biol. Chem., Feb 1996; 271: 2357.
Gretchen A. Murphy, Mol. Biol. Cell, Dec 1997; 8: 2591.
Francisco J. Nicolás, J. Cell Sci., Aug 2001; 114: 3013.
Petr Kalab, Science, Mar 2002; 295: 2452.