

## Rab7

dog, recombinant, *E. coli*

Cat. No.	Amount
PR-108	50 $\mu$ g

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -80°C

### Avoid freeze / thaw cycles

#### Form

Liquid. Supplied as a 1 mg/ml solution in 25 mM HEPES, pH 7.2, 40 mM NaCl, 5 mM DTT, 2 mM MgCl<sub>2</sub> and 10 mM GDP.

#### Activity

GTP $\gamma$ S binding: >750 mmol/mol.

#### Purity

≥90% by SDS-PAGE.

#### Description

Rab7 is a 23.5 kDa small GTPase localized in late endosomes. Rab7 regulates late endocytic membrane traffic between endosomes and lysosomes. The protein contains a C-terminal CXC motif for post-translational geranylgeranylation.

#### Selected References:

Vitelli *et al.* (1997) Role of the small GTPase RAB7 in the late endocytic pathway. *J. Biol. Chem.* **272**:4391.  
Bottger *et al.* (1996) Rab4 and Rab7 define distinct nonoverlapping endosomal compartments. *J. Biol. Chem.* **271**:29191.  
Chavrier *et al.* (1990) Localization of low-molecular-weight GTP binding-proteins to exocytic and endocytic compartments. *Cell* **62**:317.

#### Associated products available from Jena Bioscience

**Non-hydrolyzable GTP-analogs** (such as GTP $\gamma$ S, GppCp, GppNHp, NPE-caged-GTP...)

**Fluorescent GTPs** (such as MANT, ANT, TNP, or GTP labeled with a dye of your choice)

**Recombinant G-protein signaling reagents** (including GAPs, GEFs, heterotrimeric G-proteins, and many more...)

For further information please visit [www.jenabioscience.com](http://www.jenabioscience.com).