

## Rna1p (RanGAP)

*S. pombe*, recombinant, *E. coli*

Cat. No.	Amount
PR-221	50 µg

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

### Avoid freeze / thaw cycles

#### Form

Liquid. Supplied in 20 mM Tris-HCl pH 7.5, and 5 mM MgCl<sub>2</sub>.

#### Molecular Weight

43.23 kDa (386 amino acids)

#### Purity

> 95% by SDS-PAGE

### Description

Rna1p is the GTPase-activating factor (GAP) of yeast Gsp1p, an ortholog of mammalian Ran. Rna1p is located in the cytoplasm throughout the cell cycle and plays a direct role in protein import into the nucleus in *S. cerevisiae* and *S. pombe*.

### Selected References:

Hillig *et al.* (1999) The crystal structure of rna1p : a new fold for GTPase-activating protein. *Mol. Cell* **3**:781.  
Corbett *et al.* (1995) Rna1p, a Ran/TC4 GTPase activity protein, is required for nuclear import. *J. Cell Biol.* **130**:1017.

### Associated products available from Jena Bioscience

**Non-hydrolyzable GTP-analogs** (such as GTPγS, 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).**