

RIN2-RA^{GST} (residues 785-879)

Ras and Rab interactor 2 - Ras-association domain

human, recombinant, *E. coli*

Cat. No.	Amount
PR-397	500 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 25 mM HEPES pH 7.5, 50 mM NaCl, 1 mM EDTA and 50% glycerol.

Purity

> 95% by SDS-PAGE

Description

RIN2, only expressed in neurons, preferentially interacts with the GTP form of Rab5 rather than the GDP form, although it enhances the guanine exchange reaction on Rab5. Moreover, gel filtration and immunoprecipitation analyses suggest that RIN2 could form a tetramer composed of anti-parallel linkage of two parallel dimers. These results suggest that RIN2 may function as an upstream activator and/or downstream effector for Rab5 in the endocytic pathway. Additionally RIN2 interacts with several Ras proteins, it binds the GTP-loaded forms of H-Ras, TC21 and M-Ras.

Selected References:

Saito *et al.* (2002). A novel binding protein composed of homophilic tetramer exhibits unique properties for the small GTPase Rab5. *J. Biol. Chem.* **277**:3412.

Associated products available from Jena Bioscience

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

Non-hydrolyzable GTP-analogs (such as GTP γ S, GppCp, GppNHp, NPE-caged-GTP...)

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

For further information please visit www.jenabioscience.com.