

RhoA^{GST}

Ras-like GTP-binding protein, GST-tagged
human, recombinant, *E. coli*

| Cat. No. | Amount |
|----------|--------|
| PR-363 | 50 µg |

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 MgCl₂ and 10 mM GDP.

Molecular weight

19.8 kDa (178 amino acids).

Purity

≥90% by SDS-PAGE.

Description

A small GTPase of the Ras superfamily. Regulates stress fiber formation in response to growth factor stimulation.

The GST-Tag facilitates the protein's application in typical GST pull-down assays.

Selected References:

Ridley (1996) Rho: Theme and variation. *Curr. Biol.* **6**:1256.
Ridley *et al.* (1992) The small GTP-binding protein Rho regulates the assembly of focal adhesions and actin stress fibers in response to growth factors. *Cell* **70**:389.

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.