

RhoA

human, recombinant, *E. coli*

Cat. No.	Amount
PR-106	50 μ g

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 25 mM HEPES, pH 7.2, 40 mM NaCl, 5 mM DTT, 2 mM MgCl₂ and 10 mM GDP.

Purity

> 90% by SDS-PAGE.

Description

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

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 GTP γ 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.