

**RhoA**

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

Cat. No.	Amount
PR-106	50 µg

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 21.7 kDa

Accession number: AF498970

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 25 mM HEPES pH 7.5, 40 mM NaCl, 1 mM DTT, 2 mM MgCl₂ and 20 µM GDP)

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.