

## RhoC

human, recombinant, *E. coli*

Cat. No.	Amount
PR-198	50 $\mu$ 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, 2 mM DTT, 2 mM MgCl<sub>2</sub> and 10  $\mu$ M GDP.

#### Purity

> 90% by SDS-PAGE

### Description

A small GTPase of the Ras superfamily. Regulates stress fiber formation in response to growth factor stimulation. Overexpression of this gene is associated with tumor cell proliferation and metastasis.

### 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.