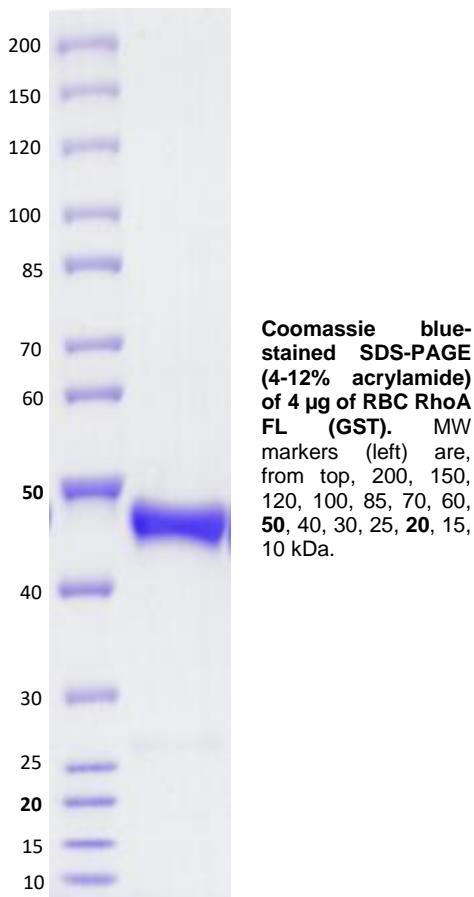


**RhoA FL (GST)****CATALOG NO.:** MSC-11-688**LOT NO.:****DESCRIPTION:** Human recombinant RhoA FL (residues 2-193; Genbank Accession # NM\_001664.4; MW = 48.5 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.**PURITY:** > 90% by SDS-PAGE.**SUPPLIED AS:** \_\_\_ µg/µL in 25 mM Tris, pH 8.0, 500 mM NaCl, 5% glycerol, 5mM MgCl<sub>2</sub>, 2 mM DTT as determined by OD<sub>595</sub>.**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice/ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

**Reaction Biology**

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com