

RhoA FL (His)

CATALOG NO.: MSC-11-687

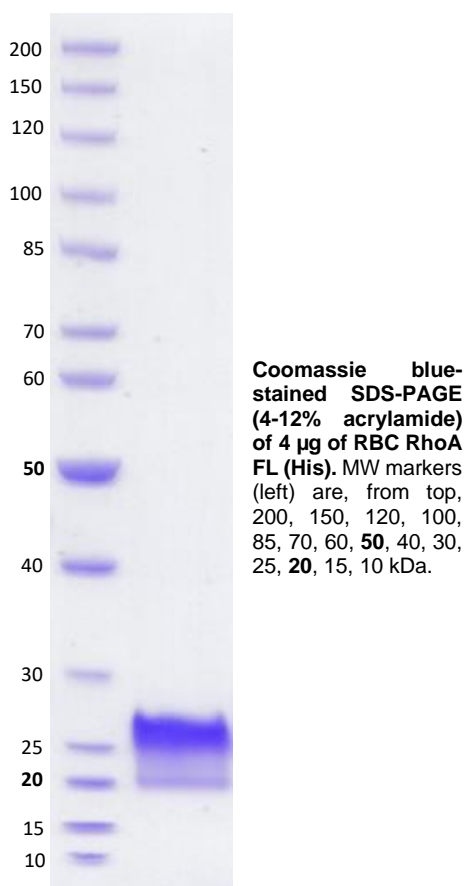
LOT NO.:

DESCRIPTION: Human recombinant RhoA FL (residues 2-193; Genbank Accession # NM_001664.4; MW = 23.8 kDa) expressed as an N-terminal His-fusion protein in *E. coli*.

PURITY: > 85% by SDS-PAGE.

SUPPLIED AS: ___ µg/µL in 25 mM Tris, pH 8.0, 500 mM NaCl, 5% glycerol, 5mM MgCl₂, 2 mM DTT as determined by OD₅₉₅.

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