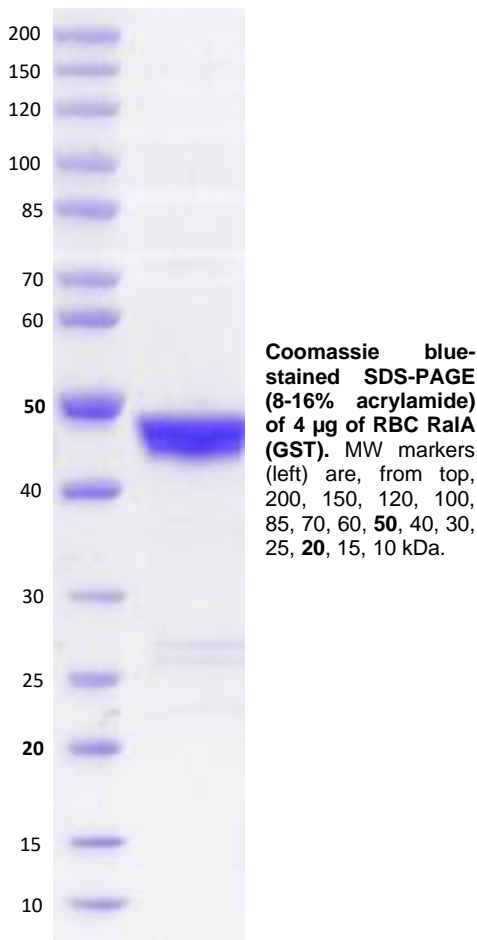


RalA FL (GST)
CATALOG NO.: MSC-11-689

LOT NO.:
DESCRIPTION: Human recombinant RalA (residues 2-206; Genbank Accession # NM_005402.4; MW = 50.3 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

PURITY: > 85% by SDS-PAGE.

SUPPLIED AS: ___ µg/µL in 50 mM Tris, pH 8.0, 300 mM NaCl, 10% glycerol, 5mM MgCl₂, 0.25 mM TCEP 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

