

HRAS (wildtype)

CATALOG NO.: MSC-11-549

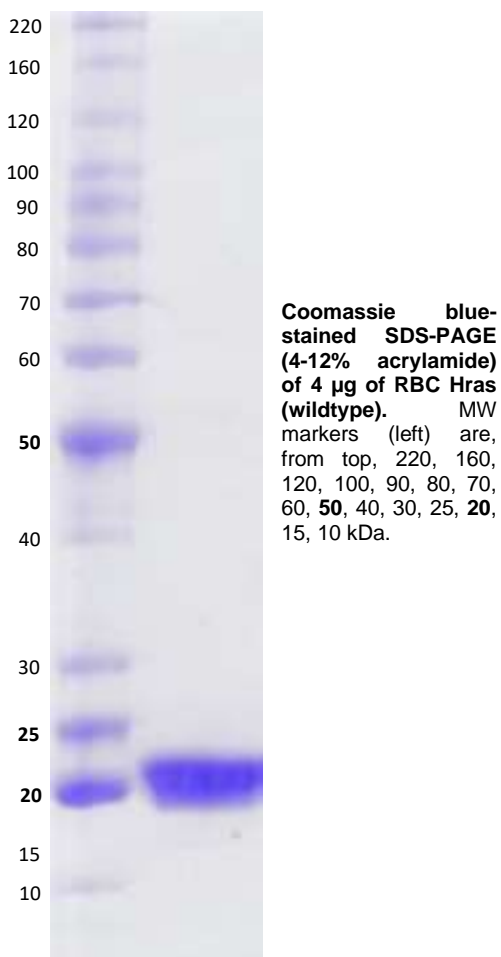
LOT NO.:

DESCRIPTION: Human recombinant HRAS (wildtype) (residues 2-169; Genbank Accession # NP_005334.1; MW = 21.3 kDa) expressed as an N-terminal His-tag protein in *E. coli*.

PURITY: > 85% by SDS-PAGE.

SUPPLIED AS: µg/µL in 25 mM Tris, pH 8.0, 100 mM NaCl, 5% glycerol, 1mM MgCl₂, 1 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