

HRAS (wildtype) (GST)

CATALOG NO.: MSC-11-551

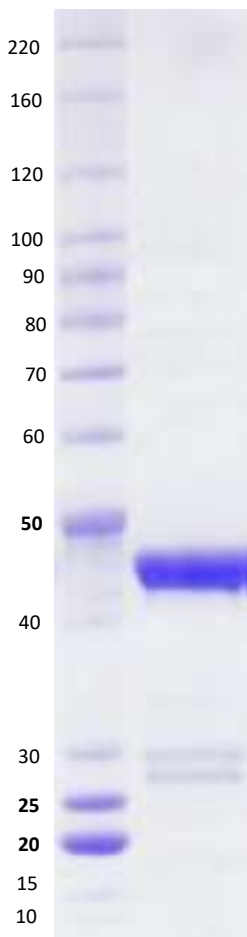
LOT NO.:

DESCRIPTION: Human recombinant HRAS (wildtype) (residues 2-169; Genbank Accession # NP_005334.1; MW = 45.99 kDa) expressed as an N-terminal GST-fusion 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.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 µg of RBC Hras (wildtype) (GST). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.

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