

**NRAS (Q61H) (GST)**

CATALOG NO.: MSC-11-605

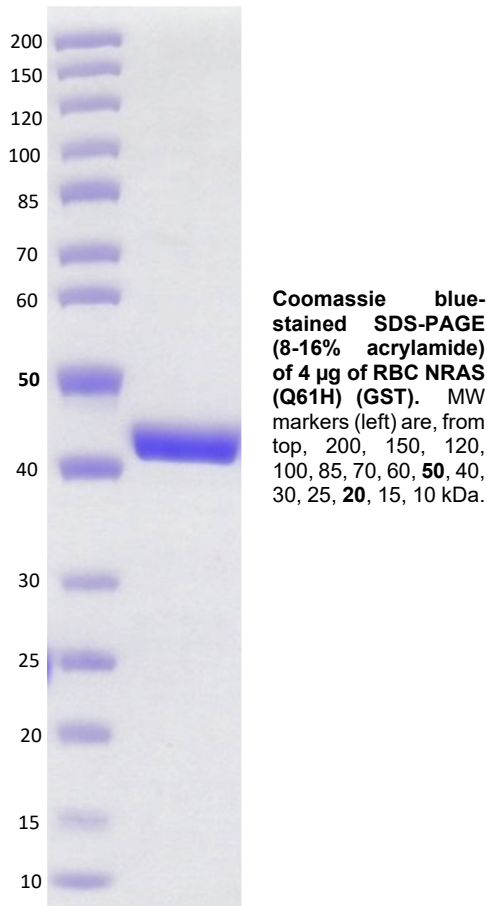
LOT NO.:

**DESCRIPTION:** Mutant human recombinant NRAS (Q61H) with histidine (H) substituted for glutamine-61 (Q61) (otherwise contains wildtype residues 2-169; Genbank Accession #NP\_002515.1; MW = 45.9 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

**PURITY:** > 95% by SDS-PAGE.

**SUPPLIED AS:** \_\_\_ µg/µL in 25 mM Tris, pH 8.0, 500 mM NaCl, 5% glycerol, 5mM MgCl<sub>2</sub>, 1 mM TCEP 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