

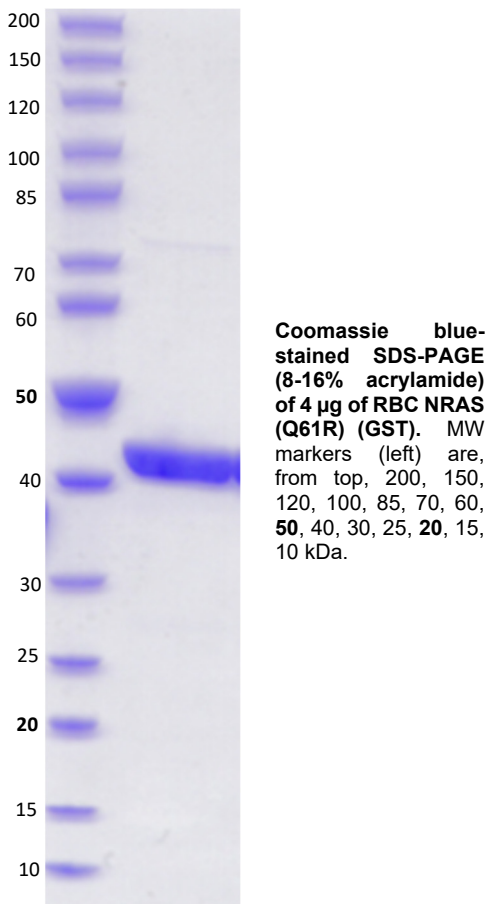
**NRAS (Q61R) (GST)****CATALOG NO.:** MSC-11-608**LOT NO.:**

**DESCRIPTION:** Mutant human recombinant NRAS (Q61R) with arginine (R) substituted for glutamine-61 (Q61) (otherwise contains wildtype residues 2-169; Genbank Accession # NP\_002515.1; MW = 46.0 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

**PURITY:** > 90% by SDS-PAGE.

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