

KRAS (Q61H)

CATALOG NO.: MSC-11-546

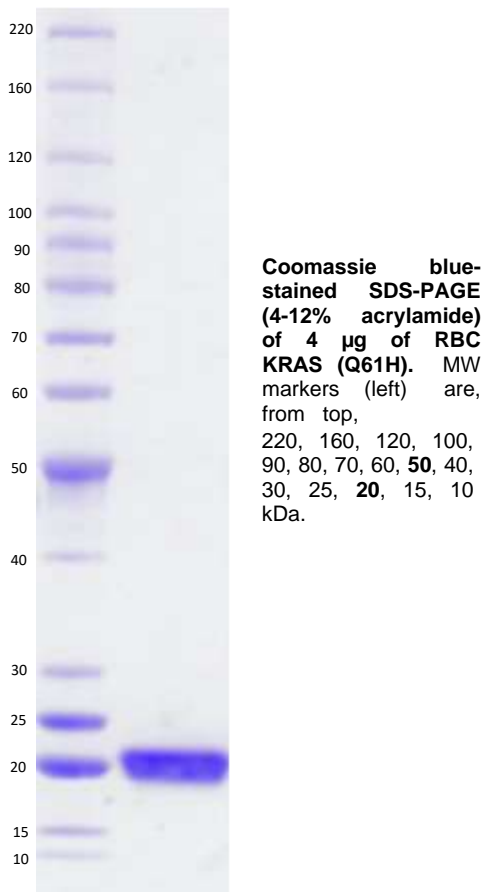
LOT NO.:

DESCRIPTION: Mutant human recombinant KRAS (Q61H) with histidine (H) substituted for glutamine-61 (Q61) (otherwise contains wildtype residues 2-169; Genbank Accession #NP_004976.2; MW = 21.3 kDa) expressed as an N-terminal His-fusion protein in *E. coli*.

PURITY: >90 % by SDS-PAGE.

SUPPLIED AS: $\mu\text{g}/\mu\text{L}$ in 25 mM Tris, pH 7.5, 150 mM NaCl, 5 mM MgCl_2 , 0.5 mM TCEP as determined by OD_{595} .

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