

KRAS (G12V) (GST)

CATALOG NO.: MSC-11-540

LOT NO.:

DESCRIPTION: Mutant human recombinant KRAS (G12V) with valine (V) substituted for glycine-12 (G12) (otherwise contains wildtype residues 2-169; Genbank Accession #NP_004976.2; MW = 46.0 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

PURITY: >85% by SDS-PAGE.

SUPPLIED AS: $\mu\text{g}/\mu\text{L}$ in 25 mM Tris, pH 8.0, 100 mM NaCl, 5% glycerol, 1 mM TCEP, 5 mM MgCl_2 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 ($<5 \mu\text{l}$) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 μg of RBC KRAS (G12V) (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.