

Biotinylated Hras (G12C)

CATALOG NO.: MSC-11-708

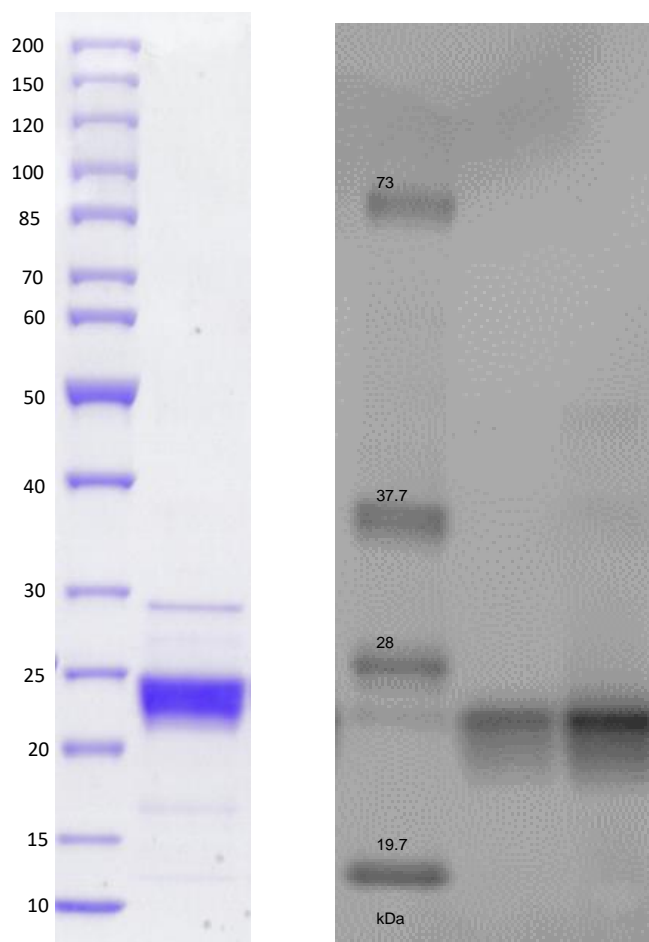
LOT NO.:

DESCRIPTION: Mutant human recombinant Hras (G12C) with cysteine (C) substituted for glycine-12 (G12) (otherwise contains wildtype residues 2-169; Genbank Accession #NP_005334.1; MW = 23.8 kDa) expressed as an N-terminal 8xHis-fusion and C-terminal BSP-tag protein in *E. coli*.

PURITY: > 85% by SDS-PAGE.

SUPPLIED AS: ___ µg/µL in 25 mM Tris-HCl pH 8.0, 500 mM NaCl, 5% glycerol, 5 mM MgCl₂, 2 mM DTT 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 (8-16% acrylamide) of 4 µg of RBC biotinylated Hras (G12C). MW markers of ladder are 200, 150, 120, 100, 85, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.

Western blot of 0.25 and 0.75 µg biotinylated Hras (G12C). Biotinylated protein was detected using IRDye 680TL conjugated streptavidin (Licor IRDye690LT 926-68031) and imaged using LI-COR Odyssey Fx. Mw markers are as indicated.

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

