

Biotinylated NRAS (wildtype)

CATALOG NO.: MSC-11-543

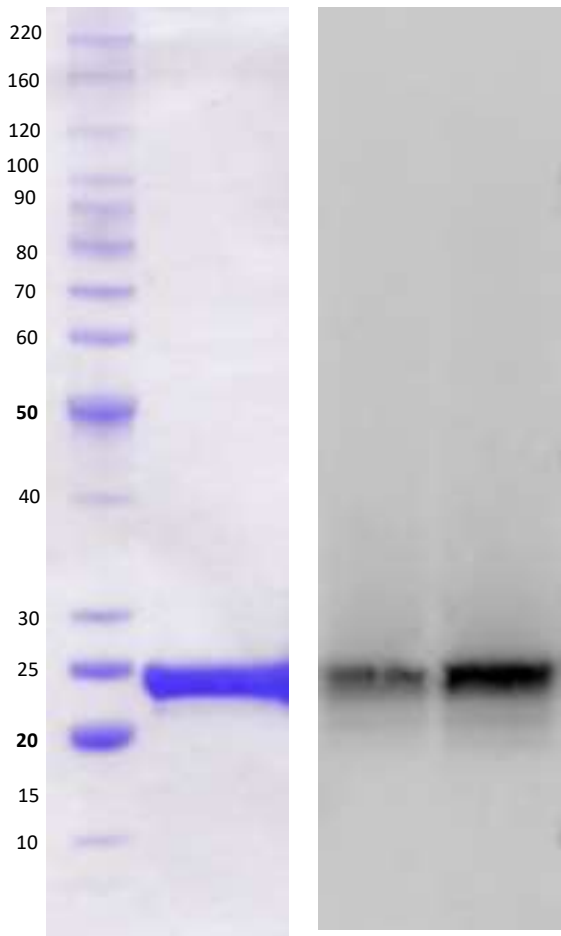
LOT NO.:

DESCRIPTION: Human recombinant NRAS (wildtype) (residues 2-169; Genbank Accession #NP_002515.1; MW = 23.7 kDa) expressed as an N-terminal 8xHis-fusion and C-terminal BSP-tag protein in *E. coli*.

PURITY: >90% by SDS-PAGE

SUPPLIED AS: µg/µL in 25 mM Tris, pH 8.0, 100 mM NaCl, 1 mM TCEP, 1 mM MgCl₂, 5% glycerol 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 (4-12% acrylamide) of 4 µg of RBC biotinylated NRAS wildtype. MW markers of ladder are indicated.

Western blot of 0.25 and 0.75µg biotinylated NRAS wildtype. Biotinylated protein was detected using IRDye 680TL conjugated streptavidin (Licor IRDye690LT 926-68031) and imaged using LI-COR Odyssey Fx.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

