

cRAF-[RBD] (GST)

CATALOG NO.: MSC-11-572

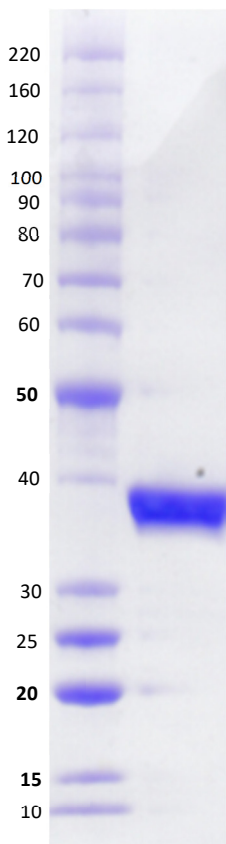
LOT NO.:

DESCRIPTION: Human recombinant cRAF-[RBD] (residues 50-140; Genbank Accession #NM_002880.3; MW = 37.18 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

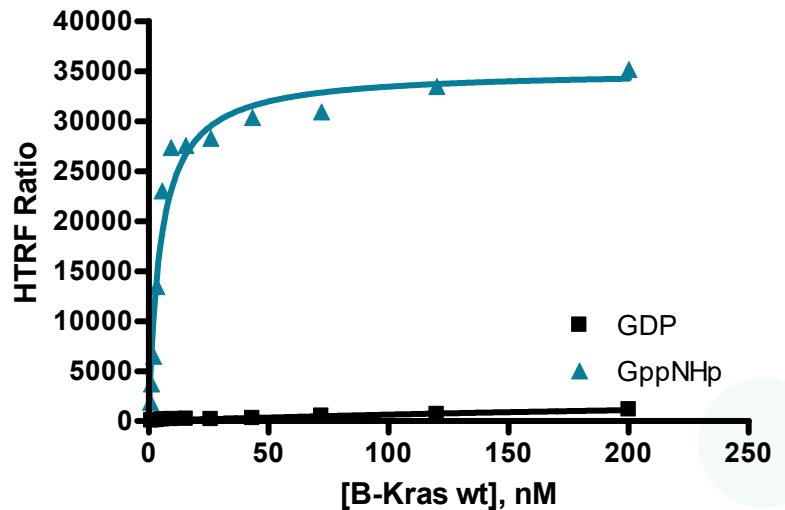
PURITY: >90% by SDS-PAGE

SUPPLIED AS: _____ µg/µL in 50 mM Tris-HCl, pH 7.5, 300 mM NaCl, 10% glycerol, 0.25 mM TCEP as determined by OD₂₈₀.

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 cRAF-[RBD] (GST). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



KRAS/cRAF PPI assay. cRAF-[RBD] binding to KRAS::GppNHp and KRAS::GDP was assessed using Perkin Elmer HTRF detection (cat #61GSTTLB and 610SAXLB). The 15µL reaction contained 5nM cRAF-[RBD], variable concentration of KRAS and detection mix. Fluorescence emission (665 and 620nm) was read using PHERAstar reader (BMGLabtech) following 2h incubation.

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