

**bRAF-[RBD] (GST)**

**CATALOG NO.:** MSC-11-574

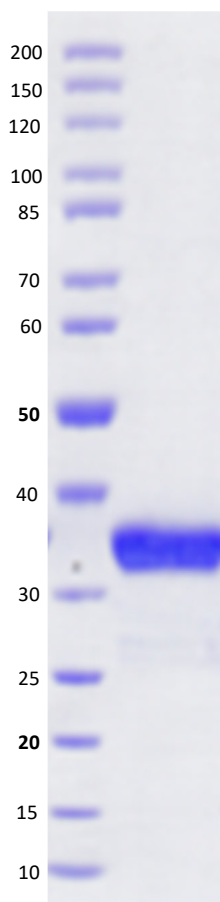
**LOT NO.:**

**DESCRIPTION:** Human recombinant bRAF-[RBD] (residues 150-237; Genbank Accession #NP\_004324.2; MW = 37.87 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

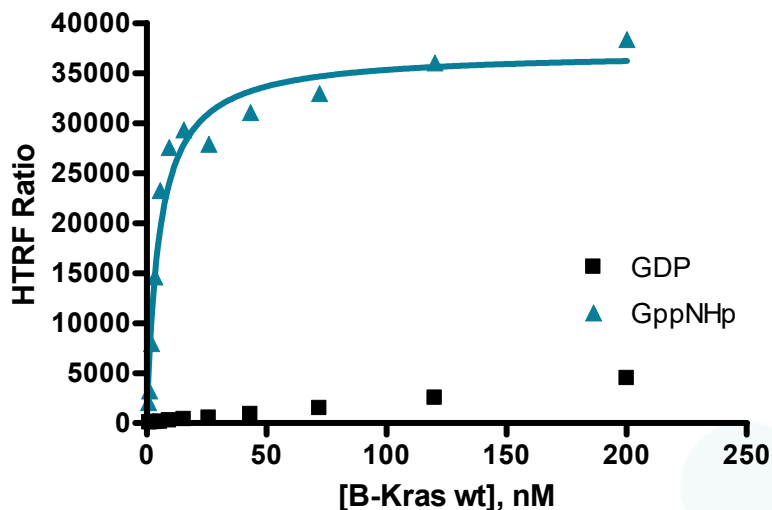
**PURITY:** >90% by SDS-PAGE

**SUPPLIED AS:** \_\_\_ µg/µL in 50 mM Tris-HCl, pH 7.5, 500 mM NaCl, 10% glycerol, 0.25 mM TCEP as determined by OD<sub>280</sub>.

**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 bRAF-[RBD] (GST). MW markers (left) are, from top, 200, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**KRAS/bRAF PPI assay.** bRAF-[RBD] binding to KRAS::GppNHp and KRAS::GDP was assessed using Perkin Elmer HTRF detection (cat #61GSTTLB and 610SAXLB). The 15µL reaction contained 5nM bRAF-[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