

Recombinant HRAS (1-186) protein

Catalog No: 81481, 81581

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 1 µg/µl

Source: Human

Buffer Contents: Recombinant HRAS (1-186) protein is supplied in 25 mM Tris 8.0, 300mM NaCl, 20% glycerol, 0.5 mM TCEP.

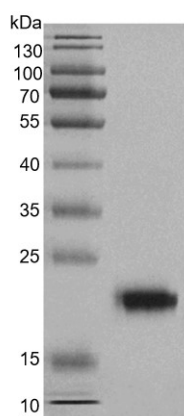
Background: **HRAS (GTPase HRas)** protein encoded by HRAS is a GTPase. **HRas** is a small G protein that belongs to the small GTPase superfamily. When HRas is combined with guanosine triphosphate, it will bind to Raf kinases such as c-Raf and further activate MAPK/ERK path. Mutations in this gene have been associated with a variety of cancers, including bladder cancer, follicular thyroid cancer, and oral squamous cell carcinoma.

Protein Details: Recombinant HRAS (1-186) protein that includes amino acids 1-186 of human HRAS protein (accession number NP_005334.1) was expressed in *E. coli* cells with an C-terminal 6xHis tag. The molecular weight of the protein is 22.06 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

HRAS (1-186)



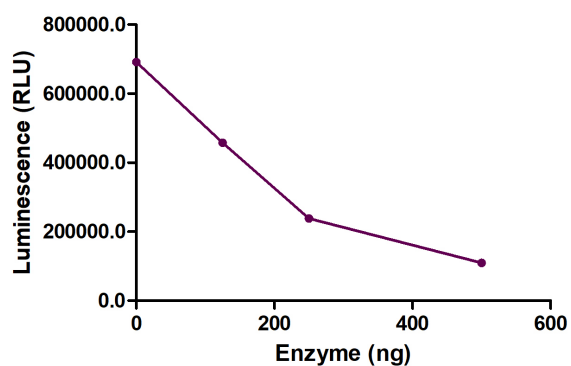
Recombinant MRM1-MBD protein

12.5% SDS-PAGE gel stained with Coomassie Blue.

MW: 22.06 kDa

Purity: >95%

HRAS (1-186) Titration



GTPase-Glo assay for HRAS (1-186) activity

1 μ M rGTP and 0.5 mM DTT was incubated with different concentrations of HRAS (1-186) in GTPase/GAP Buffer for 60 minutes at room temperature, the total reaction volume was 5 μ l. To the completed GTPase reactions, we added 5 μ l of reconstituted GTPase-Glo™ Reagent to each well and incubated the plate for 30 minutes. Then detection Reagent (10 μ l) was added, reactions were incubated for 10 minutes and luminescence was measured.