

Recombinant EEF1G protein

Catalog No: 81222, 81922

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 0.6 µg/µl

Source: Human

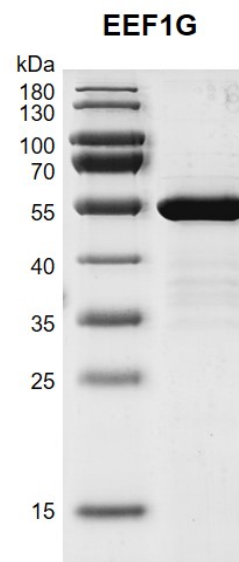
Buffer Contents: Recombinant EEF1G protein is supplied in 25 mM Tris-HCl pH 8.0, 500 mM NaCl, 5% glycerol.

Background: EEF1G (Eukaryotic Translation Elongation Factor 1 Gamma), also called EF1G or GIG35, is a subunit of the elongation factor-1 complex, which is responsible for the enzymatic delivery of aminoacyl tRNAs to the ribosome. This subunit contains an N-terminal glutathione transferase domain, which may be involved in regulating the assembly of multisubunit complexes containing this elongation factor and aminoacyl-tRNA synthetases. EEF1G probably plays a role in anchoring the complex to other cellular components.

Protein Details: Recombinant EEF1G protein was expressed in *E. coli* as the full length protein (accession number NP_001395.1) with an N-terminal 6×His tag. The molecular weight of the protein is 53.7 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 this product is shown.

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.



Recombinant EEF1G protein gel
12.5% SDS-PAGE gel
MW: 53.7 kDa
Purity: >92%