

Recombinant EEF1B2 protein

Catalog No: 81221, 81921

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 0.5 µg/µl

Source: Human

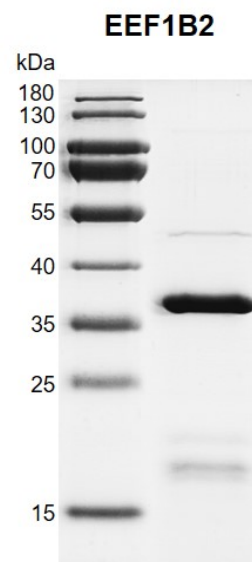
Buffer Contents: Recombinant EEF1B2 protein is supplied in 25 mM Tris-HCl pH 8.0, 500 mM NaCl, 5% glycerol.

Background: EEF1B2 (Eukaryotic Translation Elongation Factor 1 Beta 2), also called EEF1B, is specifically transcribed in fetal brain and in skeletal muscle. It is a guanine nucleotide exchange factor involved in the transfer of aminoacylated tRNAs to the ribosome. EF-1-beta (EEF1B2) and EF-1-delta (EEF1D) stimulate the exchange of GDP bound to EF-1-alpha (EEF1A1/A2) to GTP. It may also mediate accelerated synthesis of free thiol-containing proteins in response to HC-induced oxidative stress by homocysteine (HC).

Protein Details: Recombinant EEF1B2 protein was expressed in *E. coli* as the full length protein (accession number NP_001950.1) with an N-terminal 6×His-Tag. The molecular weight of the protein is 28.3 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 EEF1B2 protein gel
12.5% SDS-PAGE gel
MW: 28.3 kDa
Purity: >90%