

## Recombinant EEF1A2 protein

**Catalog No:** 81220, 81920

**Expressed In:** *E. coli*

**Quantity:** 20, 1000 µg

**Concentration:** 0.2 µg/µl

**Source:** Human

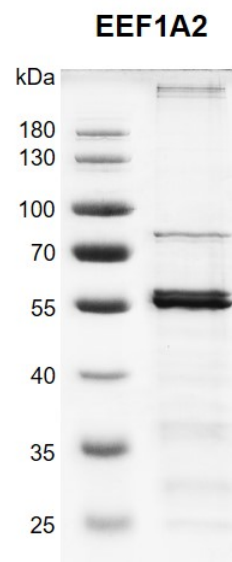
**Buffer Contents:** Recombinant EEF1A2 protein is supplied in 25 mM Tris-HCl pH 7.4, 500 mM NaCl, 10% glycerol, and 0.5 mM TCEP.

**Background:** EEF1A2 (Eukaryotic Translation Elongation Factor 1 Alpha 2), also called as EF1A or Statin, is an isoform of the alpha subunit of the elongation factor-1 complex. This complex is responsible for the enzymatic delivery of aminoacyl tRNAs to the ribosome. There are two isoforms for subunit alpha, EEF1A1 and EEF1A2. EEF1A2 promotes the GTP-dependent binding of aminoacyl-tRNA to the A-site of ribosomes during protein biosynthesis. This isoform (alpha 2) is expressed in brain, heart and skeletal muscle, and the other isoform (alpha 1) is expressed in brain, placenta, lung, liver, kidney, and pancreas. This gene may be critical in the development of ovarian cancer.

**Protein Details:** Recombinant EEF1A2 protein was expressed in *E. coli* as the full length protein (accession number NP\_001949.1) with an N-terminal 6×His tag. The molecular weight of the protein is 54 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 EEF1A2 protein gel**  
10% SDS-PAGE gel  
MW: 54 kDa  
Purity: >80%