

Recombinant eIF2 α Protein

Catalog No: 31108

Source: Human

Expressed In: Sf9 cells

Contents: 25 μ g of Recombinant eIF2 α Protein

Protein Details

Recombinant human active eIF2 α (full length) was produced in Sf9 cells and purified via sequential chromatography to >95%. Recombinant eIF2 α is phosphorylatable in vitro using either immunoprecipitated PKR or extracts from IFN γ stimulated HEK 293 cells.

Dilution of Protein

Briefly centrifuge the vial prior to opening to bring contents to the bottom of the tube. Reconstitute in PBS to a concentration of 0.2-1.0 mg/ml. Following reconstitution, supplement the protein solution with 1 mM β mercaptoethanol (final concentration) and 10% glycerol (final concentration) to prevent precipitation and store in aliquots at -80°C.

Application

Recombinant eIF2 α is suitable for Western blotting assays. The molecular weight of eIF2 α is ~36 kDa.

NOTE: The presence of Poly [d(I-C)] in buffers may affect protein functionality and should be avoided.

Storage

Lyophilized proteins can be stored at 4°C, preferably desiccated. 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 guaranteed for 6 months from date of arrival.

Research use

This product is for research use only and is not for use in diagnostic procedures.