

Recombinant I κ B α Protein

Catalog No: 31309

Source: Human

Expressed In: *E. coli*

Contents: 50 μ g of Recombinant I κ B α Protein

Protein Details

Recombinant human active I κ B α (amino acids 1-317) is produced in *E. coli* with an amino terminal polyhistidine tag and purified via sequential chromatography to >95%. Recombinant I κ B α is phosphorylatable in vitro using either recombinant active IKK α or IKK β or IKK immunoprecipitated from stimulated 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 I κ B α is suitable for Western blotting and ELISAs. The molecular weight of I κ B 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.