

Recombinant CK2 β protein

Catalog No: 31151

Source: Human

Quantity: 10 μ g

Expressed In: *E. coli*

Buffer: Protein is supplied lyophilized.

Contents: Recombinant CK2 β protein

Protein Details:

Recombinant human CK2 β was expressed in *E. coli* as a non-fusion protein. Accession number X16937. Purification by several chromatography steps to > 98% by SDS-PAGE. The beta subunit is the non-catalytic subunit of serine/threonine protein kinase CK2 and possible interacts with other protein kinases as well. The beta subunit shows no kinase activity, confers stability to the CK2 α subunit, and is involved in activity and substrate specificity.

Dilution of Protein:

The CK2 β protein is supplied lyophilized. Reconstitute in distilled water or 50% glycerol. Once reconstituted, aliquot into small fractions to avoid Freeze/Thaw cycles.

Application:

The molecular weight of the protein is 24.9 kDa.

Storage:

Some proteins are supplied lyophilized and some are supplied in solution. 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.

Guarantee:

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.