

Recombinant CK2α1 protein

Catalog No: 31149

Lot No: 01806001

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 0.4 µg/µl

Source: Human

Buffer Contents: 10 µg of Recombinant CK2α1 protein in 25 mM Tris-HCl, 500 mM NaCl, 1 mM DTT, 500 µM PMSF, and 50% glycerol, pH 8.5. Protein is supplied at 0.4 µg/µl.

Protein Details: Recombinant human CK2α1 was expressed in *E. coli* as a non-fusion protein. The accession number is M55265. The alpha subunit is the catalytic subunit of serine/threonine protein kinase CK2 and is constitutively active. This protein suitable for labeling CK2α substrates. This protein has shown to be extremely salt sensitive and highest activity is obtained without salt. Purification by several chromatography steps to > 99% by SDS-PAGE.

Application Notes: Recombinant CK2α1 is suitable for kinase assays. The molecular weight of the protein is ~45 kDa. The activity of the protein is ~ 1,320,000 units/mg with 1 unit defined as the amount of enzyme that will catalyze the transfer of 1 pmol phosphate to synthetic peptide substrate RRRDDDSDDD per minute at 37°C.

Recommended Kinase Reaction Buffer: 20 mM Tris-HCL, 25 mM β-glycerophosphate, 5 mM EGTA, and 1 mM DTT, pH 7.5.

Kinase activity may vary depending on the substrate and reaction conditions used.

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 guaranteed for 6 months from date of receipt.

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