

Recombinant PKCα protein

Catalog No: 31159

Lot No: 01806001

Expressed In: Baculovirus

Quantity: 10 µg

Concentration: 0.4 µg/µl

Source: Human

Buffer Contents: 10 µg of Recombinant PKCα protein in 20 mM NaH₂PO₄, 500 mM NaCl, 1 mM sodium orthovanadate, 5 mM NaF, 40 mM beta-glycerophosphate, 5% glycerol, pH 7.8.

Protein Details: Recombinant human PKCα protein was expressed in Sf9 cells with a C-terminal His-Tag and is suitable for labeling PKC substrates. The accession number is X52479. Purified by Ni-agarose chromatography to > 95% by SDS-PAGE.

Application Notes: Recombinant PKCα is suitable for kinase assays. The molecular weight of the protein is ~81.912 kDa. The activity of the protein is ~ 355,000 units/mg with 1 unit defined as the amount of enzyme that will catalyze the transfer of 1 pmol phosphate to PKC peptide (QKRPSQRSKYL) per minute at 30°C.

Recommended Kinase Reaction Buffer: 20 mM Tris-HCL, 25 mM β-glycerophosphate, 5 mM EGTA, 1 mM sodium orthovanadate 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.