

## Recombinant PKA protein

---

**Catalog No:** 31158

**Lot No:** 01806001

**Expressed In:** *E. coli*

**Quantity:** 50 µg

**Concentration:** 1.7 µg/µl

**Source:** Human

**Buffer Contents:** 50 µg of Recombinant PKA protein in 30 mM K-phosphate, 150 mM KCl, 1 mM EDTA, 1 mM DTT, 50% glycerol, pH 7.4.

**Protein Details:** Recombinant human PKA protein was expressed in *E. coli* with an N-terminal His-Tag and is suitable for labeling PKA substrates. The accession number is NM\_002730. Purified by Ni-agarose chromatography to > 95% by SDS-PAGE.

**Application Notes:** Recombinant PKA is suitable for kinase assays. The molecular weight of the protein is ~42.301 kDa. The activity of the protein is ~ 1,600,000 units/mg with 1 unit defined as the amount of enzyme that will catalyze the transfer of 1 pmol phosphate to PKA peptide (LRRASLG) 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.