

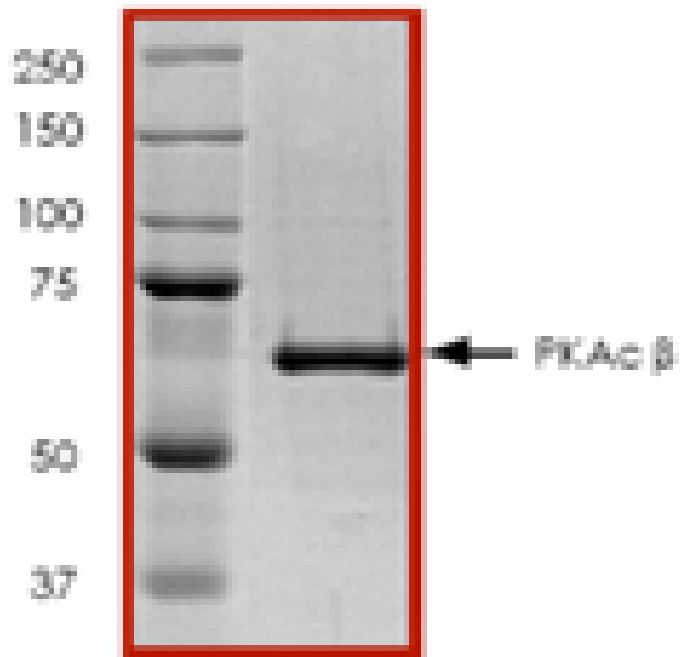
PKA Catalytic β (human), (recombinant) (GST- tag)

PKA belongs to the cAMP dependent protein kinase (PKA) family. PKA isoforms type I and II include respective dimeric R subunits RI and RII which each produce two subunits designated RI- α , RI- β , RII- α , and RII- β . The catalytic subunit also consists of more than one gene product.

Researchers cloned mammalian catalytic subunit C- α , C- β , and C- γ cDNAs. The catalytic subunit C- β belongs to the Ser/Thr protein kinase family. Activated by cAMP, PKA C- β resides in the cytoplasm (inactive holoenzyme and monomeric catalytic subunit), and translocates into the nucleus (monomeric catalytic subunit). A number of inactive tetrameric holoenzymes result from the combination of homo- or heterodimers of the different regulatory subunits associated with the two catalytic subunits. cAMP causes the dissociation of the inactive holoenzyme into a dimer of regulatory subunits bound to four cAMP and two free monomeric catalytic subunits. PKA C- β acts as a p75 neurotrophin receptor (NTR)-interacting protein, which phosphorylates p75 (NTR) at Ser304.

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis of PKA Catalytic β (human),
(recombinant).

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	cAMP-dependent protein kinase catalytic subunit β , Protein kinase A C- β
Formulation	Liquid. In 50mM Tris-HCl, pH 7.5, containing 150mM sodium chloride, 0.25mM DTT, 0.1mM EGTA, 0.1mM EDTA, 0.1mM PMSF, and 25% glycerol.
MW	~65kDa
Purity	≥80% (SDS-PAGE)
Purity Detail	Purified by SDS-PAGE / densitometry.
Source	Produced in Sf9 insect cells using an N-terminal GST-tag. Produced in a baculovirus expression system.
UniProt ID	P22694

Last modified: October 9, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)