## PKA C-α (human), (recombinant) (GSTtag)

Most of the effects of cAMP in the eukaryotic cell are mediated through the phosphorylation of target proteins on serine or threonine by the cAMP-dependent protein kinase. The inactive cAMP-dependent protein kinase is a tetramer composed of 2 regulatory and 2 catalytic subunits. The cooperative binding of 4 molecules of cAMP dissociates the enzyme in a regulatory subunit dimer and 2 free active catalytic subunits. Null mutations in PKAca leads to early postnatal death in the majority of C- $\alpha$  knock out mice.

## **Ordering Information**

Order Online »

BML-SE410-0005

5µg

Manuals, SDS & CofA

**View Online »** 

## **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 α,

Protein kinase A C-α

MW 69 kDa

Source Produced in insect cells. Full length active enzyme fused

at the N-terminus to a GST-tag. Produced in a baculovirus

expression system.

UniProt ID P17612

Last modified: May 29, 2024

