PKCα (human), (recombinant) (GSTtag)

PKC α is a member of the PKC family of serine- and threonine-specific protein kinases that can be activated by calcium and diacylglycerol. It has been reported to play roles in many different cellular processes, such as cell adhesion, cell transformation, cell cycle checkpoint, and cell volume control. Knockout studies in mice suggest that this kinase may be a fundamental regulator of cardiac contractility and Ca²⁺ handling in myocytes.

Citations: 2

View Online »

Ordering Information

Order Online »

BML-SE494-0005

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 Protein kinase Cα

Application Notes Useful for kinetic and functional studies, phosphorylation of

target substrates, drug screening.

MW 103 kDa

Source Produced in insect cells. Full length active PKCa (protein

kinase C α) is fused at the N-terminus to a GST-tag.

Produced in a baculovirus expression system.

UniProt ID P17252

Last modified: May 29, 2024



eu@enzolifesciences.com