PKCθ (human), (recombinant)

PKC θ is a member of the PKC family of serine and threonine protein kinases which can be activated by calcium and diacylglycerol.

Ordering Information

Order Online »

BML-SE517-0005

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 C θ

Application Notes Useful for kinetic and functional studies, phosphorylation of

target substrates, drug screening.

MW 110 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 Q04759

Last modified: May 29, 2024



eu@enzolifesciences.com