

PKCv (human), (recombinant) (GST- tag)

Closely related to PKC μ , PKCv has been reported to be involved in the regulation of protein trafficking in the trans-Golgi network and in the insulin-independent basal glucose uptake in L6 skeletal muscle cells.

Ordering Information

[Order Online »](#)

BML-SE349-0005	5 μ g
----------------	-----------

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 α , Protein kinase D3, PKD3

Application Notes Useful for kinetic and functional studies, phosphorylation of target substrates, drug screening.

MW 120 kDa

Source Produced in insect cells. Active, full length protein kinase C α (PKC α , protein kinase D3, PKD3) is fused at the N-terminus to a GST-tag. Produced in a baculovirus expression system.

Technical Info / Product Notes Replacement for **ADI-PPK-457**

UniProt ID O94806



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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