

# PKC $\nu$ (human), (recombinant) (GST- tag)

Closely related to PKC $\mu$ , PKC $\nu$  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.

**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 v, 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 nu (PKCv, 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

**Last modified: May 29, 2024**



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)