

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

PKC $\mu$  is a novel member of the PKC family and contains a pleckstrin homology (PH) domain. PKC $\mu$  phosphorylates HDAC7, causing it to be exported out of the nucleus and relieving Nur77 repression.

Citations: 1

[View Online »](#)

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  $\mu$ , Protein kinase D, PKD

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

**MW** 131 kDa

**Source** Produced in Sf9 insect cells. Active, full length protein kinase C mu (PKC $\mu$ , protein kinase D, PKD) is fused at the N-terminus to a GST-tag. Produced in a baculovirus expression system.

**Technical Info / Product Notes** Replacement for **ADI-PPK-450**

**UniProt ID** Q15139

**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)