

PKC α (human), (recombinant) (GST- tag)

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	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 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



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)