

PKC Lipid Activator, 10x

Catalog # L51-39

Lot # V2598-6

Product Description

10x stock solution used for assaying protein kinase activity of PKC family members. Vortex or sonicate prior to use.

Formulation

Frozen solution of buffer containing: 0.5mg/ml phosphatidylserine, 50ug/ml 1-stearoyl-2-linoleoyl-sn-glycerol, 50ug/ml 1-oleoyl-2-acetyl-sn-glycerol, 0.15% Triton X-100, 1mM DTT, 2mM CaCl₂, 20mM MOPS, pH7.2

Sterility

Buffer passed through a 0.2 μ filter membrane and packaged aseptically.

Storage and Stability

Store product at -20°C. For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Related Products

| Product Name | Catalog Number |
|----------------|----------------|
| PKC α | P61-18G |
| PKC β I | P62-10G |
| PKC β II | P63-10G |
| PKC δ | P64-10G |
| PKC ϵ | P65-10G |
| PKC γ | P66-10G |
| PKC η | P67-10G |
| PKC ι | P68-10G |
| PKC ζ | P75-10G |

PKC Lipid Activator, 10x

| | |
|--------------------|---|
| Catalog # | L51-39 |
| Lot # | V2598-6 |
| Stability | 1yr at -20°C from date of shipment |
| Storage & Shipping | Store product at -20°C. For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped frozen. |

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.