

PKC Lipid Activator, 10x

Catalog # L51-39

Lot # F733-6

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 0.2µ 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 Number L51-39

Specific Lot Number F733-6

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.