Phosphatidic acid, dipalmitoyl

PKC activator

A mitogenic lipid phosphate produced *via* the action of phospholipase D. Mediates agonist-dependent Raf-1 translocation and activation of the MAP kinase cascade.

Citations: 4

View Online »

Ordering Information

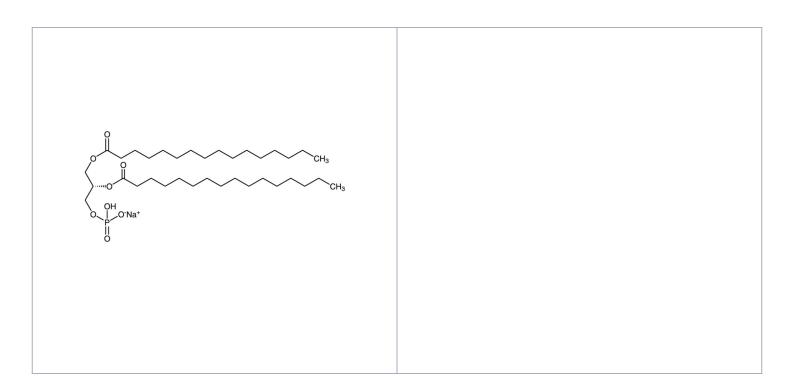
Order Online »

BML-LP103-0100

100mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up

to 3 months at -20°C.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1,2-Dipalmitoyl-sn-glycero-3-phosphate . sodium salt,

DPPA

Appearance White solid.

CAS 169051-60-9

Couple Target PKC

Couple Type Activator

Formula $C_{35}H_{68}O_8P\cdot Na$

MW 670.9

Purity ≥98% (TLC)

Solubility Fine suspension in ethanol (5mg/ml) with 15 minutes of

sonication at room temperature.

Last modified: May 29, 2024

