

# 1-Stearoyl-2-arachidonoyl-sn-glycerol

## PKC activator

Putative endogenous activator of protein kinase C produced *via* receptor-mediated inositol phospholipid hydrolysis.

Citations: 2

[View Online »](#)

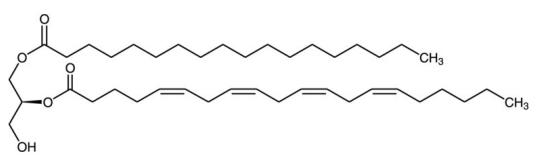
## Ordering Information

[Order Online »](#)

BML-DG130-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 2 months. Avoid contact with acids or bases. Compound may isomerize over time.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Oil or waxy solid.
CAS	65914-84-3
Couple Target	PKC
Couple Type	Activator
Formula	$C_{41}H_{72}O_5$
MW	645.0
Purity	≥95% (TLC)
Solubility	Soluble in DMSO or 100% ethanol.



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)