

# 1-O-Hexadecyl-2-O-acetyl-sn-glycerol

**PKC inhibitor**

Inhibits protein kinase C activation by diacylglycerols. Stimulates HL-60 cell differentiation. Biosynthetic precursor of PAF via remodeling and retroconversion/de novo routes. Induces delayed aggregation of rabbit platelets. Mitogenic for smooth muscle cells.

Citations: 9

[View Online »](#)

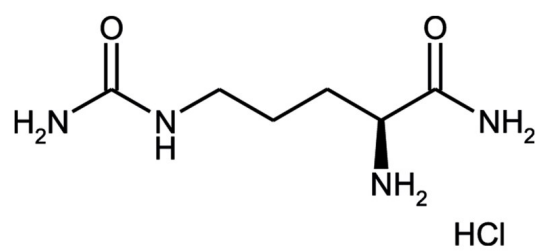
## Ordering Information

[Order Online »](#)

BML-EI150-0005	5mg
----------------	-----

## 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 at -20°C. Stock solutions are stable for up to 6 months at -20°C.
Handling	Hygroscopic. Protect from moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Waxy solid.
CAS	77133-35-8
Couple Target	PKC
Couple Type	Inhibitor
Formula	$C_{21}H_{42}O_4$
MW	358.6
Purity	≥95% (TLC)
Solubility	Soluble in 100% ethanol or acetone.

**Technical Info / Product Notes** Avoid acidic or basic conditions. Note: Product is not sterile



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)