

# 1-Oleoyl-2-acetyl-sn-glycerol

**PKC activator**

*In vitro* potent activator of protein kinase C (PKC). Stimulator of secretion and aggregation in blood platelets. Stimulates formation of actin nucleation sites. Stimulation of pp60c-src phosphorylation in EBV transformed B-cells, 470µM.

Citations: 12

[View Online »](#)

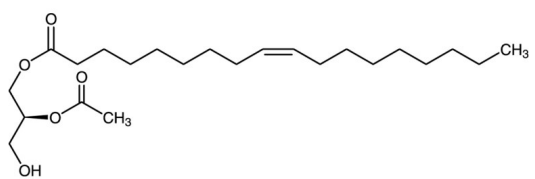
## Ordering Information

[Order Online »](#)

BML-DG125-0010	10mg
BML-DG125-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	OAG
<b>Appearance</b>	Colorless oil or waxy solid.
<b>CAS</b>	86390-77-4
<b>Couple Target</b>	PKC
<b>Couple Type</b>	Activator
<b>Formula</b>	$C_{23}H_{42}O_5$
<b>MW</b>	398.6
<b>Purity</b>	≥95% (TLC)
<b>Solubility</b>	Soluble in DMSO or ethanol.
<b>Source</b>	Synthetic.



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)