

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 »](#)

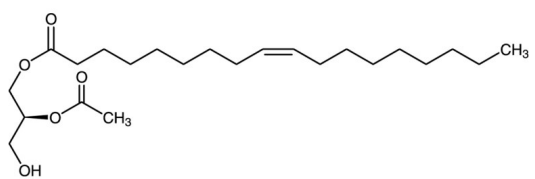
Ordering Information

[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

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

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)