

1,2-Dioctanoyl-sn-glycerol (8:0)

PKC activator

Potent activator of protein kinase C (PKC). Membrane permeable but non-physiologic DAG. Induces a discrete but transient translocation of protein kinase C.

Citations: 10

[View Online »](#)

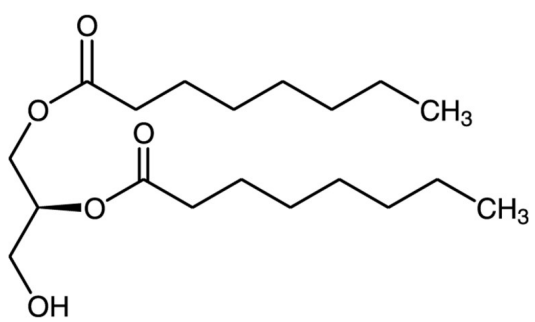
Ordering Information

[Order Online »](#)

BML-DG112-0020	20mg
-----------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	At room temperature DOG is an oil. Due to its instability, keep it cool and dry.
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DOG
Appearance	Clear colorless oil or waxy solid.
CAS	60514-48-9
Couple Target	PKC
Couple Type	Activator
Formula	$C_{19}H_{36}O_5$
Identity	Determined by 1H -NMR.
MW	344.5
Purity	≥96% (TLC)
Solubility	Soluble in DMSO or 100% ethanol.
Technical Info / Product Notes	Avoid acidic or basic conditions.

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-
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)