

7,7-

Dimethyleicosadienoic acid

Phospholipase inhibitor

Non-metabolizable analog of arachidonic or eicosadienoic acid. Cell permeable. DEDA inhibits P388D₁ cell phospholipase A₂ (PLA₂) (IC₅₀=16 μM) and snake venom PLA₂ (IC₅₀=14 μM). It inhibits human recombinant sPLA₂ (IC₅₀=5.8 μM) selectively over cPLA₂. It is also a weak inhibitor of 5-lipoxygenase (IC₅₀=56 μM, RBL-I cells). Inhibits induction of the anti-viral state in human fibroblasts induced by interferon.

Citations: 5

[View Online »](#)

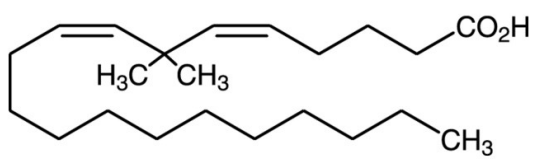
Ordering Information

[Order Online »](#)

BML-EI112-0010	10mg
BML-EI112-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at 4°C for up to 1 year. Store solutions at -20°C for up to 6 months.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DEDA
Appearance	Viscous colorless or pale yellow oil.
CAS	89560-01-0
Couple Target	Phospholipase
Couple Type	Inhibitor
Formula	$C_{22}H_{40}O_2$
MW	336.6
Purity	≥99%
Solubility	Soluble in 100% ethanol or methanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com