

Mead acid

Immunomodulatory and anti-inflammatory effects

Accumulates in tissues of essential fatty acid deficient animals. Is not converted to prostaglandins and LTB₄ but is converted to cysteinyl leukotrienes. Immunomodulatory and anti-inflammatory effects.

Citations: 3

[View Online »](#)

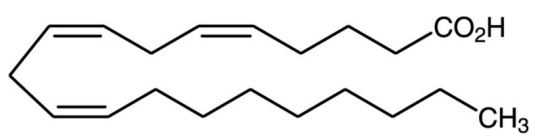
Ordering Information

[Order Online »](#)

BML-FA004-0001	1mg
BML-FA004-0010	10mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions under an inert atmosphere at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Colorless oil.
CAS	20590-32-3
Formula	$C_{20}H_{34}O_2$
MW	306.5
Purity	≥97% (HPLC)
Solubility	Soluble in ethanol, DMSO, acetone, ethyl acetate or hexane.

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)