

# Mead acid

## Immunomodulatory and anti-inflammatory effects

Accumulates in tissues of essential fatty acid deficient animals. Is not converted to prostaglandins and  $\text{LTB}_4$  but is converted to cysteinyl leukotrienes. Immunomodulatory and anti-inflammatory effects.

Citations: 3

[View Online »](#)

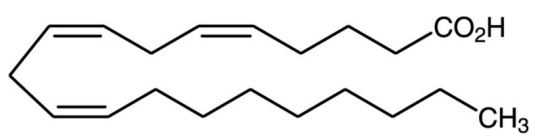
## Ordering Information

[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	Clear solution.
CAS	20590-32-3
Formula	$C_{20}H_{34}O_2$
Formulation	10 mg/mL EtOH.
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-  
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)