

# 17-Octadecynoic acid

Fatty acid  $\omega$ -hydroxylase inhibitor

Potent suicide inhibitor of LTB<sub>4</sub>  $\omega$ -oxidase (IC<sub>50</sub> < 5  $\mu$ M). Potent and selective inhibitor of fatty acid  $\omega$ -hydroxylase (IC<sub>50</sub> < 100 nM). Blocks biosynthesis of 20-HETE.

Citations: 3

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-EI103-0010

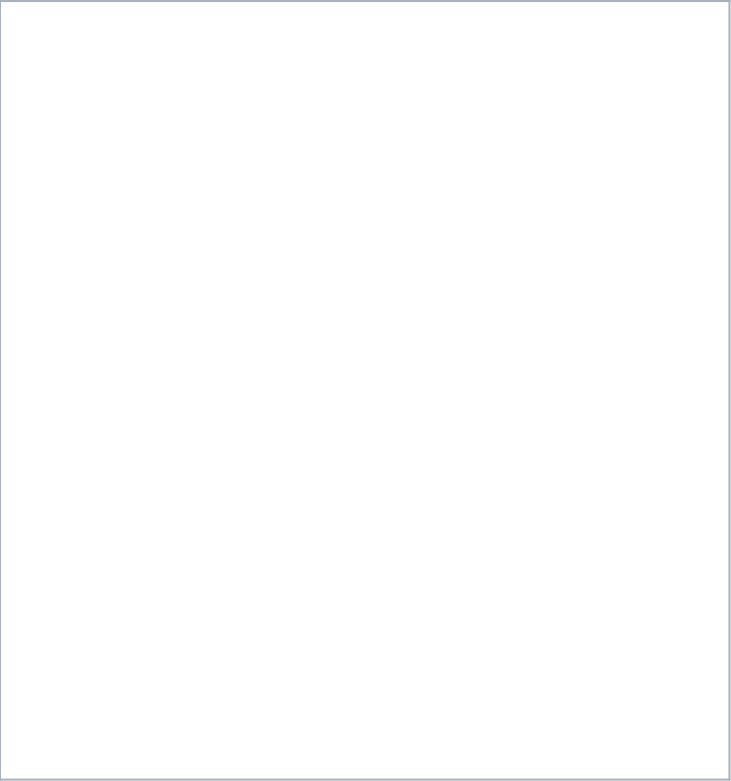
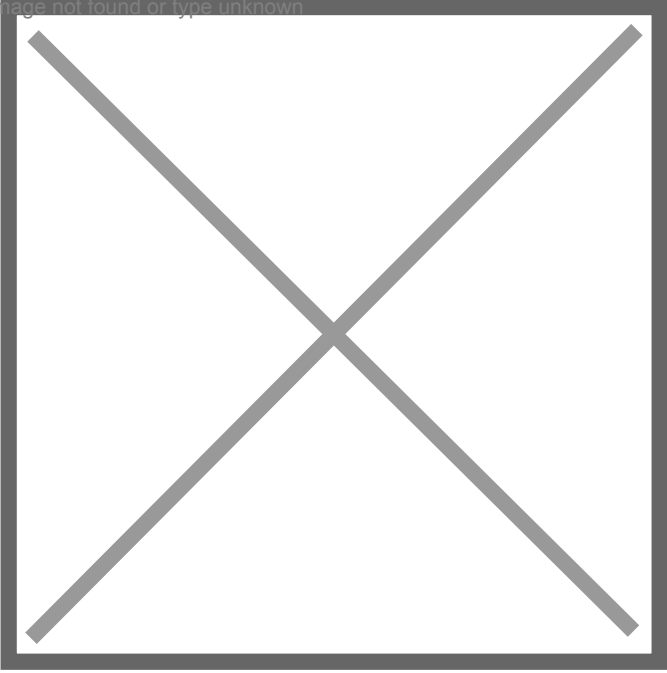
10mg

Manuals, SDS & CofA

[View Online »](#)



Image not found or type unknown



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	White solid.
CAS	34450-18-5
Couple Target	Fatty acid omega-hydroxylase, Leukotriene oxidase
Couple Type	Inhibitor
Formula	$C_{18}H_{32}O_2$
MW	280.5
MeltingPoint	67-68°C
Purity	≥98% (HPLC)
Solubility	Soluble in ethanol (25mg/ml) or DMSO (25mg/ml).

Last modified: May 29, 2024

Image not found or type not known  
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)