

(±)9-Hydroxyoctadeca- 10E,12Z-dienoic acid

Ultra-pure

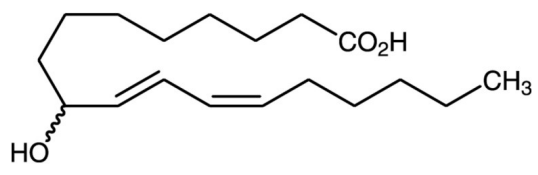
Ordering Information

[Order Online »](#)

BML-HO309-0050	50µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-80°C
Shipping	Dry Ice – Dangerous Good

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	(±)9-HODE
Appearance	Clear to light-yellow liquid.
CAS	98524-19-7
Formula	$C_{18}H_{32}O_3$
Formulation	Solution in ethanol.
MW	296.4
Purity	≥99% (HPLC)
Source	Synthetic.

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-be@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