

13(S)- Hydroxyoctadeca- 9Z,11E-dienoic acid

Inhibits thromboxane production in human platelets

Biosynthesis in rabbit peritoneal serous membranes, atherosclerotic plaques, healthy blood vessels, psoriatic skin, endothelial cells. Stimulates prostacyclin production by endothelial cells. Inhibits thromboxane production in human platelets. Inhibits platelet and tumor cell adhesion to endothelium. Antagonizes 12-HETE-induced tumor cell adhesion to endothelium. Stimulates DNA synthesis in BALB/c 3T3 fibroblasts.

Citations: 7

[View Online »](#)

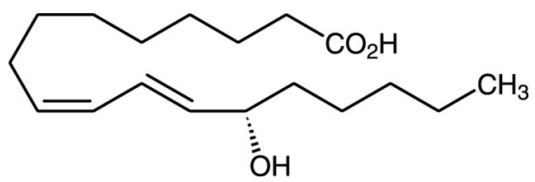
Ordering Information

[Order Online »](#)

BML-HO013-0050	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

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 13(S)-HODE

Appearance Clear to light-yellow liquid.

CAS 29623-28-7

Formula $C_{18}H_{32}O_3$

Formulation Solution in ethanol.

MW 296.5

Purity ≥99% (HPLC)

Source Synthetic.

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)