

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

Inhibits prostacyclin synthase

Suppresses activation of platelets. Induces endothelial injury. Constricts isolated coronary arteries. Inhibits prostacyclin synthase.

Citations: 3

[View Online »](#)

Ordering Information

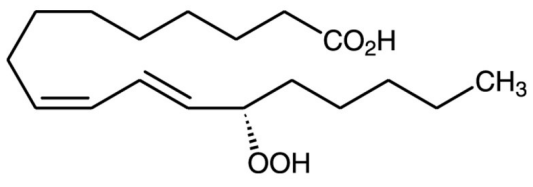
[Order Online »](#)

BML-HP013-0050

50µg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Open ampule and aliquot into suitable tubes or vials. Store at -80°C or colder. For experiments with cells or tissues, aliquots of ethanol solutions may be added directly to aqueous media. For HPLC, aliquots of ethanol solutions should be evaporated to dryness under a stream of nitrogen and HPLC mobile phase added to the desired concentration.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	13(S)-HPODE
CAS	33964-75-9
Couple Target	Prostacyclin synthase
Couple Type	Inhibitor
Formulation	Oil dissolved in ethanol.
MW	312.5
Purity	≥98%

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)