

15(S)- Hydroperoxyeicosa- 5Z,8Z,11Z,13E,17Z- pentaenoic acid

Ultra-pure
Primary metabolite of eicosapentaenoic acid *via* the 15-lipoxygenase pathway.

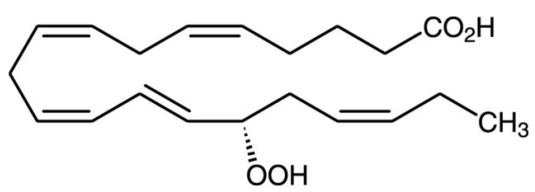
Ordering Information

[Order Online »](#)

BML-HP515-0050	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 15(S)- HPEPE

Appearance Clear to light-yellow liquid.

CAS 125992-60-1

Formula $C_{20}H_{30}O_4$

Formulation Solution in ethanol.

MW 334.5

Purity ≥98% (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)