

15(S)-Hydroxyeicosa-5Z,8Z,11Z,13E-tetraenoic acid (HETE)

Ultra-pure

Biosynthesis in human neutrophils, eosinophils and reticulocytes. Most abundant eicosanoid in human bronchi. Regulates T lymphocyte mitogenesis. Induces immunosuppression. Stimulates migration of capillary endothelial cells and neovascularization.

Citations: 8

[View Online »](#)

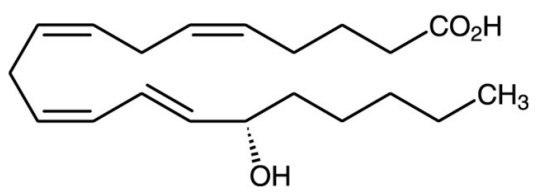
Ordering Information

[Order Online »](#)

BML-H015-1000	1mg
BML-H015-0050	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 15(S)-HETE

Appearance Liquid.

CAS 54845-95-3

Formulation Oil dissolved in ethanol.

MW 320.5

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)