

12(S)-Hydroxyeicosa-5Z,8Z,10E,14Z-tetraenoic acid (HETE)

Ultra-pure

Biosynthesis by platelet 12-lipoxygenase. Chemotactic and chemokinetic for PMNL² and for vascular smooth muscle cells. Antiproliferative effect on smooth muscle cells. Second messenger in angiotensin induced aldosterone production. Induces endothelial cell retraction and tumor cell adhesion.

Citations: 1

[View Online »](#)

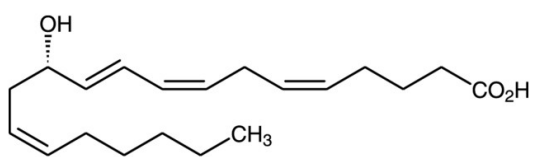
Ordering Information

[Order Online »](#)

BML-H012-0050	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 12(S)-HETE

Appearance Liquid.

CAS 54397-83-0

Formula $C_{20}H_{32}O_3$

Formulation Liquid. Solution in ethanol.

MW 320.5

Purity ≥98% (HPLC)

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)