

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

ATPase inhibitor

Biosynthesis in psoriatic skin and corneal epithelial cells via cytochrome P-450. More potent than 12(S)-HETE in inducing human leukocyte chemokinesis and aggregation. Proinflammatory effects in human skin. Inhibits renal, cardiac and corneal Na^+,K^+ -ATPase. Reduces intraocular pressure.

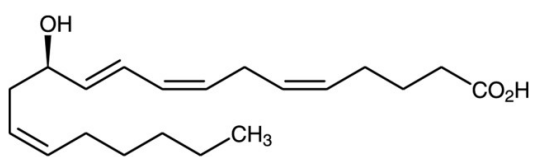
Ordering Information

[Order Online »](#)

BML-H112-0050	50µg
---------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -80°C for up to 1 year.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	12(R)-HETE
CAS	82337-46-0
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{20}H_{32}O_3$
Formulation	Liquid. Solution in ethanol.
MW	320.5
Purity	≥99%

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)