

8(S)-Hydroxyeicosa- 5Z,9E,11Z,14Z- tetraenoic acid (HETE)

Activator of PKC and PPAR α

Biosynthesis in human neutrophils, phorbol ester-treated mouse epidermis and by hepatic microsomal cytochrome P-450. Chemotactic for neutrophils. Activator of PPAR α . Activates mouse skin protein kinase C (PKC).

Citations: 4

[View Online »](#)

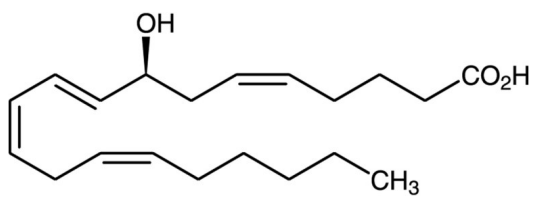
Ordering Information

[Order Online »](#)

BML-H008-0050	50 μ g
---------------	------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	8(S)-HETE
CAS	98462-03-4
Couple Target	PKC, PPAR
Couple Type	Activator
Formula	$C_{20}H_{32}O_3$
Formulation	Liquid. In ethanol.
MW	320.5
Purity	≥99% (HPLC)

Technical Info / Product Notes

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.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com