

5(S),6(R)- Dihydroxyeicosa- 7E,9E,11Z,14Z- tetraenoic acid

(DIHETE)

Synthesis via enzymatic and nonenzymatic hydrolysis of LTA₄. Induces contraction of guinea pig lung parenchymal strips. Partial agonist at LTD₄ receptors. Potent inhibitor of LTB₄ induced neutrophil migration (IC₅₀ = 8nM).

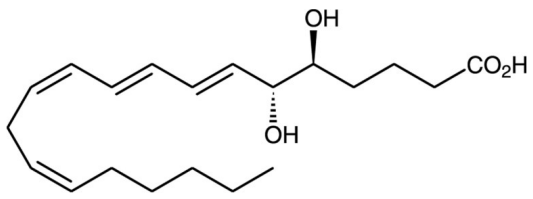
Prostanoid receptor ligand

Citations: 2

[View Online »](#)

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 5(S),6(R)-DiHETE

Appearance Colorless liquid.

CAS 82948-88-7

Couple Target Leukotriene receptor, Prostanoid receptor

Couple Type Ligand

Formula $C_{20}H_{32}O_4$

Formulation Oil dissolved in ethanol.

MW 336.5

Purity ≥99% (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)