

20-Hydroxy-leukotriene B₄

Primary LTB₄ metabolite produced by ω oxidation in PMNL retaining some chemotactic activity. ω oxidation is blocked by 17-ODYA.

Citations: 4

[View Online »](#)

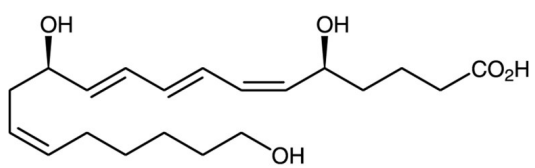
Ordering Information

[Order Online »](#)

BML-HLB004-0050	50 μ g
-----------------	------------

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 20-Hydroxy-LTB4

Appearance Clear to light-yellow liquid.

CAS 79516-82-8

Formula $C_{20}H_{32}O_5$

Formulation Solution in ethanol.

MW 352.5

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)