

Leukotriene A4 methyl ester

Substrate for LTA₄ hydrolase and LTC₄ synthase

Unstable intermediate in the biosynthesis of LTB₄ and LTC₄. The naturally occurring free acid is a substrate for LTA₄ hydrolase and LTC₄ synthase and plays a central role in transcellular leukotriene and lipoxin biosynthesis. Mobilizes Ca²⁺ in human neutrophils. Supplied as the methyl ester for greater stability.

Citations: 4

[View Online »](#)

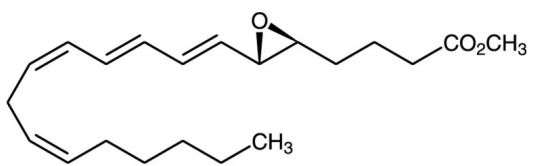
Ordering Information

[Order Online »](#)

BML-LA004-0050	50µg
----------------	------

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	LTA4-ME
Appearance	Liquid.
CAS	73466-12-3
Couple Target	Leukotriene hydrolase, Leukotriene synthase, Prostanoid receptor
Couple Type	Ligand, Substrate
Formula	$C_{21}H_{32}O_3$
Formulation	Oil dissolved in 2% triethylamine/hexane.
MW	332.5
Purity	≥95% (HPLC) (remainder 11-trans isomer)
Source	Synthetic.

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