

# Leukotriene A4 methyl ester

Substrate for LTA<sub>4</sub> hydrolase and LTC<sub>4</sub> synthase

Unstable intermediate in the biosynthesis of LTB<sub>4</sub> and LTC<sub>4</sub>. The naturally occurring free acid is a substrate for LTA<sub>4</sub> hydrolase and LTC<sub>4</sub> synthase and plays a central role in transcellular leukotriene and lipoxin biosynthesis. Mobilizes Ca<sup>2+</sup> in human neutrophils. Supplied as the methyl ester for greater stability.

Citations: 4

[View Online »](#)

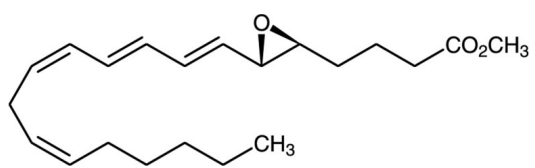
## Ordering Information

[Order Online »](#)

BML-LA004-1000	1mg
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.



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)