

N-Acetyl-leukotriene E4

Major biliary metabolite of cysteinyl leukotrienes in the rat. Approximately 100-fold less active than LTC₄. Equiactive with LTE₄ in guinea pig lung parenchyma assay.

Citations: 5

[View Online »](#)

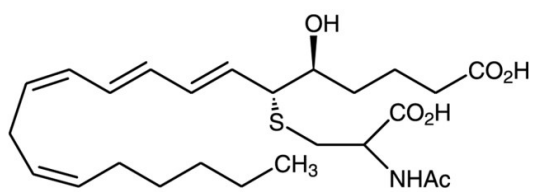
Ordering Information

[Order Online »](#)

BML-LEA004-0050	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-Acetyl-LTE4
Appearance	Liquid.
CAS	80115-95-3
Couple Target	Leukotriene receptor, Prostanoid receptor
Couple Type	Ligand
Formula	$C_{25}H_{39}NO_6S$
Formulation	Solution in ethanol.
MW	481.7
Purity	≥98% (HPLC)
Solubility	Soluble in ethanol.

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-
fr@enzolifesciences.com](mailto:info-fr@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)