

Leukotriene E4

Prostanoid receptor ligand

Less potent than LTD₄ in smooth muscle contractile activity and mucus secretion in airways. Reduces myocardial contractility and coronary blood flow. Induces long lasting hypotension and airway hyperresponsiveness. Partial agonist at LTD₄ receptors. LTE₄ solely activates a CysLT₁ receptor. Biosynthesis from LTD₄ by a dipeptidase.

Citations: 7

[View Online »](#)

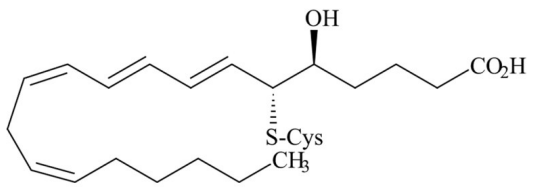
Ordering Information

[Order Online »](#)

BML-LE004-1000	1mg
BML-LE004-0050	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	LTE4
Appearance	Liquid.
CAS	75715-89-8
Couple Target	Leukotriene receptor, Prostanoid receptor
Couple Type	Ligand
Formula	$C_{23}H_{37}NO_5S$
Formulation	Solution in ethanol.
MW	439.6
Purity	≥98% (HPLC)

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