

Leukotriene C4

Prostanoid receptor ligand

A cysteinyl leukotriene which acts as potent agonist for contraction of smooth muscle. Induces microvascular permeability and mucus secretion in airways. Enhances growth of human airway epithelial cells. Reduces myocardial contractility and coronary blood flow. Induces long lasting hypotension. Stimulates angiogenesis. Stimulates LH8, LHRH9 and prolactin release. Produced in specific cell types or by transcellular biosynthesis and rapidly metabolized to LTD4 and LTE4. Transported by the multidrug resistance protein (MRP; $K_m=35\text{nM}$).

Citations: 9

[View Online »](#)

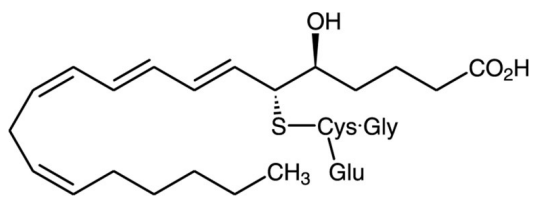
Ordering Information

[Order Online »](#)

BML-LC004-0050	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. Protect from exposure to air. Open ampule and aliquot into suitable tubes or vials. Store at -80°C or colder. For experiments with cells or tissues, aliquots of stock solutions may be evaporated to near dryness under a stream of nitrogen and dissolved in ethanol. Stock solutions may be used directly for HPLC analysis.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	LTC ₄ , 5S-Hydroxy-6R-(S-glutathionyl)-7E,9E,11Z,14Z-eicosatetraenoic acid
Appearance	Colorless solution.
CAS	72025-60-6
Couple Target	Leukotriene receptor, Prostanoid receptor
Couple Type	Ligand
Formula	C ₃₀ H ₄₇ N ₃ O ₉ S
Formulation	A solution in ethanol.
MI	14: 5455
MW	625.8
Purity	≥95% (HPLC)
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-
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)