

# Leukotriene E4

## Prostanoid receptor ligand

Less potent than LTD<sub>4</sub> 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 LTD<sub>4</sub> receptors. LTE<sub>4</sub> solely activates a CysLT<sub>1</sub> receptor. Biosynthesis from LTD<sub>4</sub> by a dipeptidase.

Citations: 7

[View Online »](#)

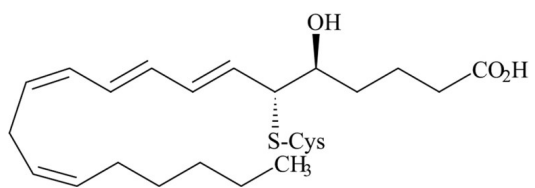
## Ordering Information

[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](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)