

# Leukotriene D4

## Prostanoid receptor ligand

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. Biosynthesis from LTC<sub>4</sub> by  $\gamma$ -glutamyltranspeptidase.

Citations: 5

[View Online »](#)

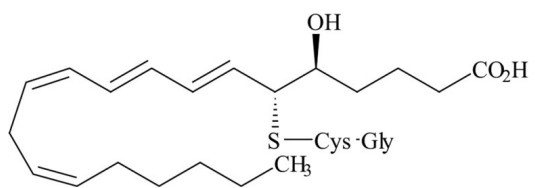
## Ordering Information

[Order Online »](#)

BML-LD004-0050	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.
Long Term Storage	-80°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	LTD4
Appearance	Liquid.
CAS	73836-78-9
Couple Target	Leukotriene receptor, Prostanoid receptor
Couple Type	Ligand
Formula	$C_{25}H_{40}N_2O_6S$
Formulation	Solution in methanol/ammonium acetate buffer, 70:30, pH 5.6.
MW	496.7
Purity	≥98% (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-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)