

Leukotriene B₄-dimethylamide

Prostanoid receptor ligand

Leukotriene B₄ antagonist. Inhibits binding of [³H]-LTB₄ to guinea pig lung membranes ($K_i=0.13\text{ }\mu\text{M}$). Antagonizes the degranulation response to LTB₄ in rabbit PMNL ($K_d=0.2\text{ }\mu\text{M}$) acting, most likely, via receptor desensitization. Partial agonist at higher concentrations (5 μM).

Citations: 1

[View Online »](#)

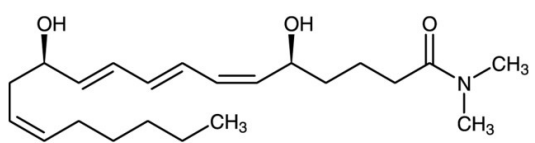
Ordering Information

[Order Online »](#)

BML-EI102-0050	50 μg
----------------	------------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store as supplied (-80°C) for up to two years.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Liquid.
CAS	83024-92-4
Couple Target	Leukotriene, Prostanoid receptor
Couple Type	Inhibitor, Ligand
Formula	$C_{22}H_{37}NO_3$
Formulation	Solution in ethanol.
MW	363.5
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-
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)