

U-75302

Prostanoid receptor ligand

Selective leukotriene B₄ antagonist, K_i=159 nM for inhibition of [³H]-LTB₄ binding to guinea pig lung membranes.

Citations: 3

[View Online »](#)

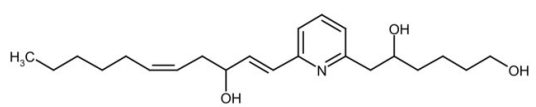
Ordering Information

[Order Online »](#)

BML-RA102-0100	100µg
----------------	-------

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. Store solutions at -80°C for up to 1 month.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Colorless oil.
CAS	119477-85-9
Couple Target	Leukotriene, Prostanoid receptor
Couple Type	Inhibitor, Ligand
Formula	$C_{22}H_{35}NO_3$
MW	361.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25 mg/ml) or ethanol (25 mg/ml).

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