

BML-111

Lipoxin receptor ligand

Potent lipoxin A₄ agonist, equiactive as LXA₄ in inhibition of LTB₄ induced PMNL chemotaxis (IC₅₀=5 nM).

Citations: 4

[View Online »](#)

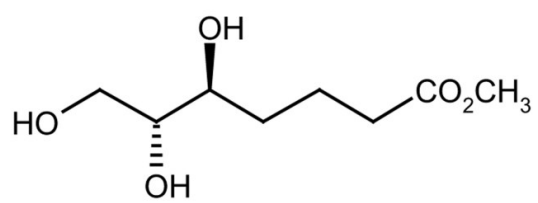
Ordering Information

[Order Online »](#)

BML-RA108-0001	1mg
BML-RA108-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5(S),6(R),7-Trihydroxyheptanoic acid methyl ester
Appearance	White solid.
Couple Target	Lipoxin receptor
Couple Type	Ligand
Formula	$C_8H_{16}O_5$
MW	192.2
Purity	≥99%
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/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