

Lipoxygenase inhibitor

Selective 12-lipoxygenase inhibitor. IC_{50} =0.063 μ M, 3.33 μ M and 1.89 μ M for 12-, 15- and 5-LO respectively in isolated enzyme preparations.

Citations: 3

[View Online »](#)

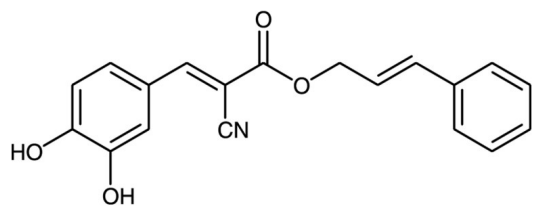
Ordering Information

[Order Online »](#)

BML-EI211-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store under nitrogen after opening. Product is subject to oxidation. Make solutions fresh daily.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

CAS 132465-11-3

Couple Target Lipoxygenase

Couple Type Inhibitor

MW 321.3

Purity ≥98%

Solubility Soluble in ethanol (5mg/ml) or DMSO (5mg/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