

Lipoxygenase inhibitor

Selective 12-lipoxygenase inhibitor. IC₅₀=0.063 μM, 3.33 μM and 1.89 μM for 12-, 15- and 5-LO respectively in isolated enzyme preparations.

Citations: 3

[View Online »](#)

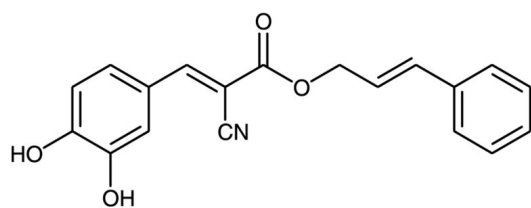
Ordering Information

[Order Online »](#)

BML-EI211-0010	10mg
BML-EI211-0050	50mg

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).



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)