

5,8,11-Eicosatriynoic acid

Lipoxygenase inhibitor

Inhibits lipoxygenases (ID_{50} =3, 20 and 20 μ M for 5-, 12- and 15-LO respectively) selectively over cyclooxygenase (ID_{50} =50 μ M) in whole cells. Also inhibits hepatic glutathione S-transferase and leukotriene C₄ biosynthesis.

Citations: 5

[View Online »](#)

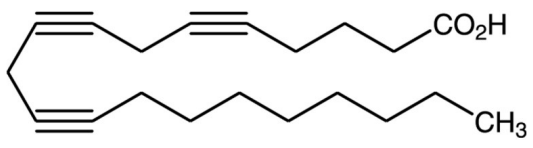
Ordering Information

[Order Online »](#)

BML-ET003-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at 0-4°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light and oxygen. Keep under inert gas.
Long Term Storage	+4°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	ETI
Appearance	Off-white solid.
CAS	13488-22-7
Couple Target	GST, Lipoxygenase
Couple Type	Inhibitor
Formula	$C_{20}H_{28}O_2$
MW	300.4
Purity	≥98% (TLC)
Solubility	Soluble in ethanol (25mg/ml), DMSO (25mg/ml) or dimethyl formamide.

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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)