

5,8,11,14- Eicosatetraynoic acid

Arachidonic acid uptake inhibitor

Inhibits arachidonic acid uptake. Inhibits cyclooxygenase ($ID_{50}=8\text{ }\mu\text{M}$) and all lipoxygenases ($ID_{50}=10, 0.3, \text{ and } 0.2\text{ }\mu\text{M}$ for 5-, 12- and 15-LO respectively) in whole cells. Inhibits PLA_2 and cytochrome P-450. Modulates Ca^{2+} entry into cells. Stimulates luteinizing hormone release from cultured pituitary cells. PPAR agonist.

Citations: 13

[View Online »](#)

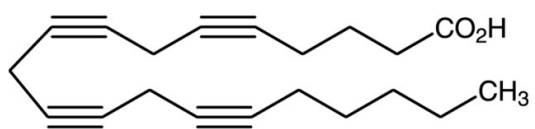
Ordering Information

[Order Online »](#)

BML-ET004-0020	20mg
BML-ET004-0100	100mg

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. Make solutions fresh daily.
Handling	Protect from light and oxygen.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ETYA
Appearance	Off-white solid.
CAS	1191-85-1
Couple Target	COX, Cytochrome P450, Lipoyxygenase, PLA, PPAR
Couple Type	Activator, Inhibitor
Formula	$C_{20}H_{24}O_2$
MW	296.5
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol, methanol, DMSO, acetone or acetonitrile.

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-
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)