

Linoleamide

Unsaturated analog of endogenous sleep inducing lipid

Unsaturated analog of endogenous sleep-inducing lipid.

Citations: 2

[View Online »](#)

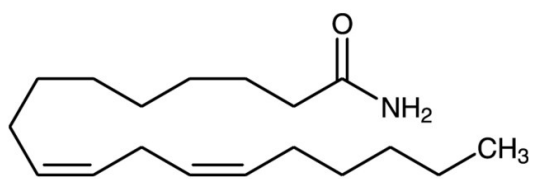
Ordering Information

[Order Online »](#)

BML-FA023-0010	10mg
BML-FA023-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from oxygen and light. After opening, protect from oxidation. Dissolve in solvent purged with nitrogen or argon. Store in tightly sealed container under nitrogen or argon at -80°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	3072-13-7
Formula	$C_{18}H_{33}NO$
MW	279.5
MeltingPoint	57-58°C
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol or in DMSO.

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-
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)