

N-Arachidonylglycine

FAAH inhibitor

Endogenous anandamide-like compound. Lacks affinity for CB₁ receptors ($K_i > 10 \mu\text{M}$), TRPV1 ($\text{EC}_{50} > 10 \mu\text{M}$) and anandamide uptake ($\text{IC}_{50} > 50 \mu\text{M}$), but inhibits fatty acid amide hydrolase (FAAH) ($\text{IC}_{50} = 8.5 \mu\text{M} - 50 \mu\text{M}$, depending on cell type and species). Displays anti-inflammatory and analgesic activity. Identified as an insulin secretagogue in a primary β -cell-based functional assay. Represents a new class of "lipoamino acids".

Citations: 6

[View Online »](#)

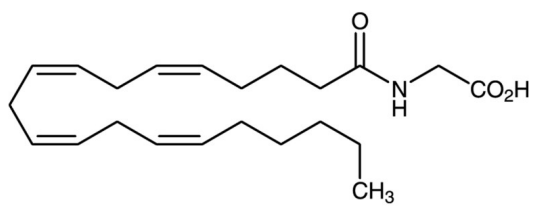
Ordering Information

[Order Online »](#)

BML-FA029-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from air.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White waxy solid.
CAS	179113-91-8
Couple Target	FAAH
Couple Type	Inhibitor
Formula	$C_{22}H_{35}NO_3$
MW	361.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO or 100% ethanol. 2mg/ml soluble in PBS (pH 7.2).
Source	Synthetic.



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)