

Arachidonylcyclopropylamide

Cannabinoid receptor ligand

High affinity selective agonist for the cannabinoid CB₁ receptor (K_i = 2.2 nM; CB₂ receptor K_i = 0.7 μ M). Inhibits forskolin-induced cAMP accumulation, increases binding of GTP γ S to cerebellar membranes, inhibits electrically-evoked contractions of mouse *vas deferens* and production of hypothermia in mice. Active *in vivo*.

Citations: 3

[View Online »](#)

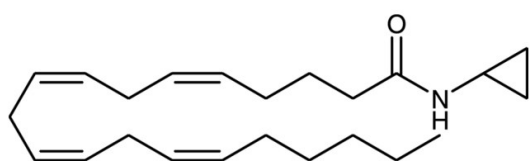
Ordering Information

[Order Online »](#)

BML-FA031-0010	10mg
BML-FA031-0050	50mg

Manuals, SDS & CoFA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Pale yellow oil.
CAS	229021-64-1
Couple Target	Cannabinoid receptor
Couple Type	Ligand
Formula	$C_{23}H_{37}NO$
MW	343.6
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml), 100% ethanol (25mg/ml) or dimethyl formamide. 1mg/ml soluble in a 1:1 solution of ethanol:PBS (pH 7.2) (dilute the ethanol solution with PBS).

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)