

Virodhamine

Cannabinoid receptor ligand

Endogenous cannabinoid which has been found to occur at concentrations 2 to 9 fold greater than anandamide in peripheral tissue. It acts as a full agonist at the CB₂ receptor and an antagonist at the CB₁ receptor.

Citations: 2

[View Online »](#)

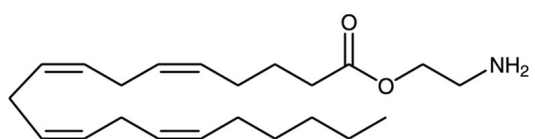
Ordering Information

[Order Online »](#)

BML-FA036-0005	5mg
----------------	-----

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 -80°C and under an inert atmosphere.
Handling	Protect from oxygen. Keep under inert gas.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	O-AEA, O-Arachidonoylethanolamine, Arachidonic acid-(2-aminoethyl)-ester, O-(2-aminoethyl)-5Z,8Z,11Z,14Z-eicosatetraenoic acid ester
Appearance	Yellow to brown oil.
CAS	443129-35-9
Couple Target	Cannabinoid receptor
Couple Type	Ligand
Formula	$C_{22}H_{37}NO_2$
MW	347.5
Purity	≥95% (TLC)
Solubility	Soluble in DMSO (45mg/ml), dimethyl formamide (40mg/ml) or ethanol (20mg/ml)

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)