

Virodhamine

Cannabinoid receptor ligand

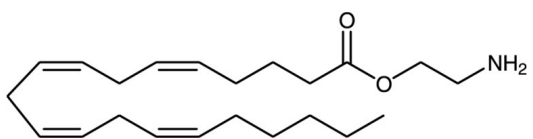
Endogenous cannabinoid which has been found to occur at concentrations 2 to 9 fold greater than anandamide in peripheral tissue. It acts as a full agonist at the CB₂ receptor and an antagonist at the CB₁ receptor.

Citations: 2

[View Online »](#)

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-Arachidonylethanolamine, Arachidonic acid-(2-aminoethyl)-ester, O-(2-aminoethyl)-5Z,8Z,11Z,14Z-icosatetraenoic 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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com