

# Anandamide

## Endogenous CB<sub>1</sub> receptor ligand

Endogenous [1] ligand for the CB<sub>1</sub> receptor (CB<sub>1</sub>: K<sub>i</sub>=52nm; CB<sub>2</sub>: K<sub>i</sub>=1930nm [2]) and TRPV1 (K<sub>i</sub>=5.78μM [3,4]). Inhibits NF-κB activation through direct binding to IKKβ [5] and induces apoptosis independently of cannabinoid or vanilloid receptors [6]. Activates the MAP kinase (MAPK/ERK) signalling pathway [7].

Citations: 30

[View Online »](#)

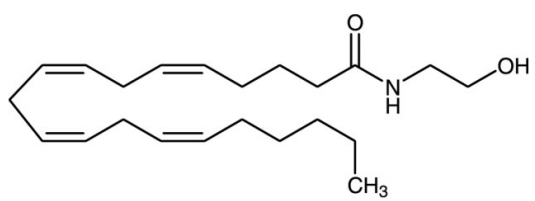
## Ordering Information

[Order Online »](#)

|                |      |
|----------------|------|
| BML-FA017-0005 | 5mg  |
| BML-FA017-0025 | 25mg |

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

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                                                                                              |
|------------------|--------------------------------------------------------------------------------------------------------------|
| Alternative Name | AEA, N-Arachidonoyl ethanolamine, (all-Z)-N-(2-Hydroxyethyl)-5,8,11,14-eicosatetraenamide                    |
| Appearance       | Colorless oil.                                                                                               |
| CAS              | 94421-68-8                                                                                                   |
| Couple Target    | Cannabinoid receptor, IKK, MAPK, TRP channel                                                                 |
| Couple Type      | Activator, Inhibitor, Ligand                                                                                 |
| Formula          | $C_{22}H_{37}NO_2$                                                                                           |
| MI               | 14: 624                                                                                                      |
| MW               | 347.54                                                                                                       |
| Purity           | ≥98% (TLC)                                                                                                   |
| Solubility       | Soluble in ethanol (e.g. at a concentration of 50 mg/ml). It is also soluble in DMSO and dimethyl formamide. |

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)