

Methanandamide

Cannabinoid receptor ligand

Amidase resistant cannabinoid receptor (CB) agonist (CB₁: K_i=20nM; CB₂: K_i=815nM). The most potent of the series of methyl-anandamides. About 4-fold higher binding affinity for cannabinoid receptor CB₁ than anandamide (Prod. No. BML-FA017) in the presence of PMSF. Does also bind to TRPV1 (K_i=4.67μM). Inhibits motor behavior resembling the effects of THC rather than anandamide.

Citations: 9

[View Online »](#)

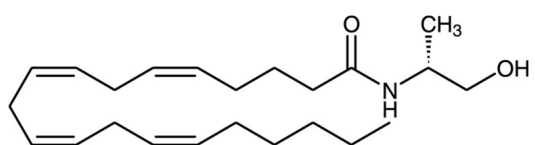
Ordering Information

[Order Online »](#)

BML-FA021-0005	5mg
BML-FA021-0025	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, at -80°C. Stock solutions are stable for up to 1 month at -20°C.
Handling	Protect from light and oxygen. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(R)-(+)-Arachidonyl-1'-hydroxy-2'-propylamide, AM 356, N-(2-Hydroxy-1R-methylethyl)-5Z,8Z,11Z,14Z-eicosatetraenamide
Appearance	Colorless oil.
CAS	157182-49-5
Couple Target	Cannabinoid receptor
Couple Type	Ligand
Formula	$C_{23}H_{39}NO_2$
MW	361.6
Purity	≥98% (TLC)
Solubility	Soluble in ethanol or DMSO (10mg/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