

2-

Arachidonoylglycerol

Cannabinoid receptor ligand

Endogenous ligand for the CB₁ receptor (CB₁: K_i=58.3nM); CB₂: K_i>3μM) present in high levels in the CNS.

Citations: 16

[View Online »](#)

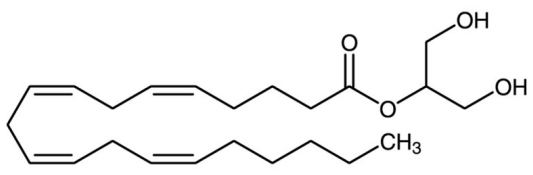
Ordering Information

[Order Online »](#)

BML-FA028-0005	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 -70°C. Stock solutions are stable for up to 3 months when stored at -70°C under an atmosphere of nitrogen or argon.
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	2-AG
Appearance	Colorless liquid.
CAS	53847-30-6
Couple Target	Cannabinoid receptor
Couple Type	Ligand
Formula	$C_{23}H_{38}O_4$
Formulation	10 mg/mL in Acetonitrile.
Identity	Determined by IR.
MW	378.6
Purity	≥95% (TLC) (as a 9:1 mixture of the 2-AG and 1-AG).
Solubility	Soluble in Acetonitrile, DMSO or 100% ethanol.

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-
be@enzolifesciences.com](mailto:info-be@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)