

Arachidonyl-2-chloroethylamide

Cannabinoid receptor ligand

Non-pungent, potent and highly selective agonist of the cannabinoid CB₁ receptor ($K_i=1.4\text{nM}$). Displays >1'400-fold selectivity over CB₂ receptors ($K_i=3.1\mu\text{M}$). Active *in vivo*.

Citations: 4

[View Online »](#)

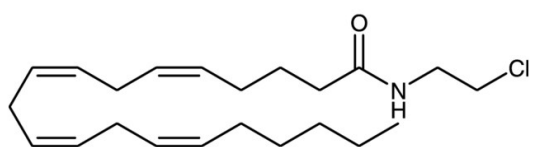
Ordering Information

[Order Online »](#)

BML-FA032-0050	50mg
----------------	------

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 -20°C. Stock solutions are stable for up to 3 months at -20°C when stored under an inert atmosphere.
Handling	Protect from air. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ACEA, N-(2-Chloroethyl)-5Z,8Z,11Z,14Z-eicosatetraenamide
Appearance	Pale yellow oil.
CAS	220556-69-4
Couple Target	Cannabinoid receptor
Couple Type	Ligand
Formula	$C_{22}H_{36}NOCl$
MW	366.0
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or ethanol (25mg/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)