

Noladin ether

Cannabinoid receptor ligand

Endogenous cannabinoid agonist isolated from porcine brain. Binds specifically to CB₁ ($K_i=21.2\text{nM}$) and CB₂ ($K_i=>3\mu\text{M}$) receptors. Causes sedation hypothermia, intestinal immobility and mild antinociception in mice. Chemically stable, but less potent than 2-arachidonoylglycerol (Prod. No. BML-FA028) with an endogenous half-life time of hours rather than minutes.

Citations: 10

[View Online »](#)

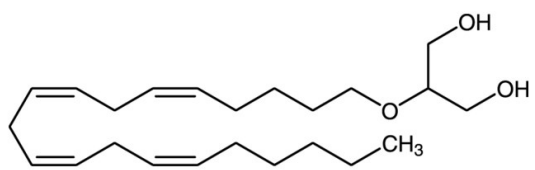
Ordering Information

[Order Online »](#)

BML-FA033-0005	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 -20°C. Stock solutions are stable for up to 3 months at -20°C when stored under an atmosphere of argon or nitrogen.
Handling	Protect from oxygen. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-Arachidonyl glyceryl ether, 2-AG ether, 5Z,8Z,11Z,14Z-Eicosatetraen-2-glyceryl ether
Appearance	Yellow oil.
CAS	222723-55-9
Couple Target	Cannabinoid receptor
Couple Type	Ligand
Formula	$C_{23}H_{40}O_3$
MW	364.6
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (40mg/ml) or ethanol (25mg/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