

Arvanil

TRPV1 channel activator

Arvanil is a non-pungent capsaicin analog with antinociceptive and anti-inflammatory effects *in vivo*. It is an agonist at TRPV1 as well as CB1, but not CB2.

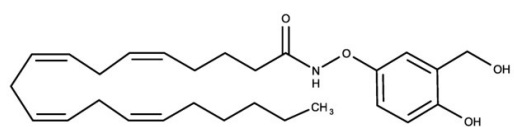
Ordering Information

[Order Online »](#)

ENZ-CHM442-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-(3-Methoxy-4-hydroxy-benzyl)-arachidonamide

Appearance Colorless or slightly yellow viscous oil.

CAS 128007-31-8

Couple Type Activator

Formula $C_{28}H_{41}NO_3$

Identity Determined by NMR.

MW 439.64

Purity ≥98% (TLC)

Solubility Soluble in DMSO (50 mg/ml).

Last modified: January 24, 2025



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)