

VDM-11

Anandamide transporter inhibitor

Selective and potent inhibitor of the anandamide membrane transporter (AMT) (neuronal (C6 glioma cells): $IC_{50}=11\mu M$; non-neuronal (rat RBL-2H3 cells): $IC_{50}=7.2\mu M$) with negligible action at TRPV1 and weak action at CB₁ receptor (CB₁: $K_i=5\mu M$; CB₂: $K_i>5\mu M$) and FAAH ($IC_{50}=3.7\mu M$ / $IC_{50}>50\mu M$ (N18TG2 cell membranes)).

Citations: 7

[View Online »](#)

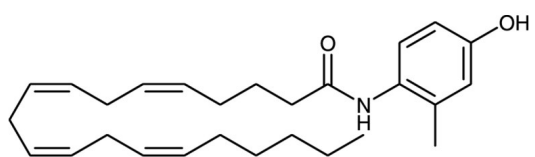
Ordering Information

[Order Online »](#)

BML-CR111-0050	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.
Handling	Keep under inert gas.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-Arachidonoyl-4-hydroxy-2-methylaniline, N-(4-Hydroxy-2-methylphenyl)arachidonoylamide
Appearance	Tan oil.
CAS	313998-81-1
Couple Target	Anandamide transporter
Couple Type	Inhibitor
Formula	$C_{27}H_{39}NO_2$
MW	409.6
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (1.5mg/ml) or ethanol (1mg/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)