

N- Arachidonoyldopamine

Activator of CB₁ receptor and TRPV1

Endogenous, specific ligand for the CB₁ receptor (CB₁: K_i=250nM; CB₂: K_i=12μM) [1] and TRPV1 [2] (EC₅₀=50nM) found in nervous tissues [3, 4]. Immunosuppressant inhibiting T cell proliferation and phosphorylation of the NF-κB p65 subunit [5]. Potent vasorelaxant [6] and inhibitor of HIV-1 [7].

Citations: 10

[View Online »](#)

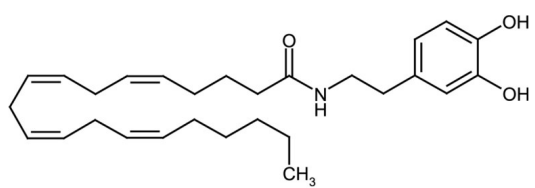
Ordering Information

[Order Online »](#)

BML-VR104-0025	25mg
----------------	------

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	Protect from oxygen. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-Arachidonoyl-3-hydroxytyramine, NADA, AA-DA
Appearance	Viscous oil.
CAS	199875-69-9
Couple Target	Cannabinoid receptor, TRP channel
Couple Type	Activator, Ligand
Formula	$C_{28}H_{41}NO_3$
MW	439.6
Purity	≥98% (TLC)
Solubility	Soluble in DMSO.

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)