

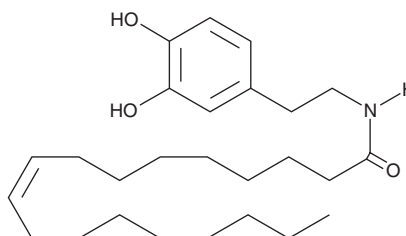
PRODUCT INFORMATION



N-Oleoyl Dopamine

Item No. 10115

CAS Registry No.: 105955-11-1
Formal Name: N-[2-(3,4-dihydroxyphenyl)ethyl]-
9Z-octadecenamide
Synonym: ODA
MF: C₂₆H₄₃NO₃
FW: 417.6
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 283 nm



Laboratory Procedures

For long term storage, we suggest that N-oleoyl dopamine (ODA) be stored as supplied at -20°C. It should be stable for at least one year.

ODA is supplied as a crystalline solid. A stock solution may be made by dissolving the ODA in the solvent of choice. ODA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of ODA in these solvents is approximately 50, 20, and 30 mg/ml, respectively.

ODA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ODA should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. ODA has a solubility of approximately 500 µg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ODA is a selective, endogenous vanilloid receptor 1 (VR₁) agonist isolated from bovine brain.¹ Structurally, it is the amide of oleic acid and dopamine and is therefore a "hybrid" analog which incorporates components of both the anadamide-like and dopamine neurotransmitter pathways. ODA binds to the human recombinant VR₁ with a K_i of 36 nM making it equipotent to capsaicin and slightly more potent than N-arachidonoyl dopamine in this assay.¹ It causes hyperalgesia and nocifensive behavior that is blocked by the VR₁ antagonist iodo-resiniferatoxin. ODA is selective for VR₁ based on observations that it has weak affinity for the rat CB₁ receptor (K_i of 1.6 µM) and is a very weak inhibitor of FAAH. ODA is also a potent inhibitor of 5-lipoxygenase from rat basophilic leukemia-1 (RBL-1) cells, with a IC₅₀ of 7.5 nM.^{2,3}

References

1. Chu, C.J., Huang, S.M., De Petrocellis, L., *et al.* N-oleoyldopamine, a novel endogenous capsaicin-like lipid that produces hyperalgesia. *J. Biol. Chem.* **278**(16), 13633-13639 (2003).
2. Tseng, C.-F., Iwakami, S., Mikajiri, A., *et al.* Inhibition of *in vitro* prostaglandin and leukotriene biosyntheses by cinnamoyl-β-phenethylamine and N-acyldopamine derivatives. *Chem. Pharm. Bull.* **40**(2), 396-400 (1992).
3. Iwakami, S., Shibuya, M., Tseng, C.-F., *et al.* Inhibition of arachidonate 5-lipoxygenase by phenolic compounds. *Chem. Pharm. Bull.* **34**, 3960-3963 (1986).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/10/2015

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM