

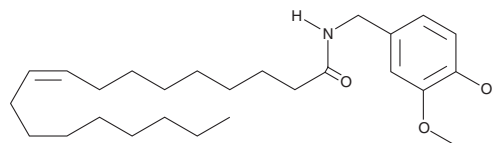
PRODUCT INFORMATION



Olvanil

Item No. 90262

CAS Registry No.: 58493-49-5
Formal Name: N-[(4-hydroxy-3-methoxyphenyl)methyl]-9Z-octadecenamide
Synonyms: NE 19550, N-Vanillyloleamide
MF: C₂₆H₄₃NO₃
FW: 417.6
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 281 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Olvanil is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of olvanil in these solvents is approximately 13 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of olvanil is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. For greater aqueous solubility, olvanil can be directly dissolved in 0.1 M Na₂CO₃ (1 mg/ml) and then diluted with PBS (pH 7.2) to achieve the desired concentration or pH. We do not recommend storing the aqueous solution for more than one day.

Description

Olvanil is a structural analog of capsaicin, which is the noxious active component of hot peppers of the *Capsicum* genus. It is the amide of vanillylamine and oleic acid. Olvanil acts as an agonist at the vanilloid receptor, VR₁, inducing desensitization analgesia in rat and mouse models of pain.¹ Olvanil has complex interactions with the cannabinoid system, in that it potentiates the agonist activity of endogenous cannabinoids by inhibiting the reuptake of arachidonyl ethanolamide. Olvanil is a more potent reuptake inhibitor than AM404, which is commonly used for this purpose (50% inhibition of reuptake at 10 μM versus 12% for AM404 at the same dose).² Olvanil is also a CB₁ agonist, but does not bind to CB₂ receptors. The IC₅₀ for FAAH is 20 μM. The overall activity of olvanil in most models is that of an analgesic, but it is unclear how these effects are mediated by VR₁, the CB₁ receptor, or other components of the endogenous pain sensation system.

References

1. Janusz, J.M., Buckwalter, B.L., Young, P.A., *et al.* Vanilloids. 1. Analogs of capsaicin with antinociceptive and antiinflammatory activity. *J. Med. Chem.* **36**, 2595-2604 (1993).
2. Di Marzo, V., Bisogno, T., Melck, D., *et al.* Interactions between synthetic vanilloids and the endogenous cannabinoid system. *FEBS Lett.* **436**, 449-454 (1998).
3. Glaser, S.T., Abumrad, N.A., Fatade, F., *et al.* Evidence against the presence of an anandamide transporter. *Proc. Natl. Acad. Sci. USA* **100**(7), 4269-4274 (2003).

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, 01/17/2023

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