

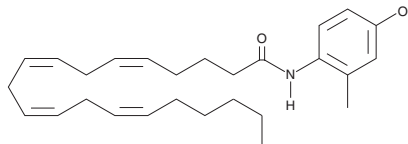
# PRODUCT INFORMATION



## VDM11

Item No. 10006731

**CAS Registry No.:** 313998-81-1  
**Formal Name:** N-(4-hydroxy-2-methylphenyl)-5Z,8Z,11Z,14Z-eicosatetraenamide  
**MF:** C<sub>27</sub>H<sub>39</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 409.6  
**Purity:** ≥98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥1 year  
**UV/Vis. λmax:** 202, 231 nm



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

### Laboratory Procedures

VDM11 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 VDM11 in these solvents is approximately 20 mg/ml.

VDM11 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of VDM11 should be diluted with the aqueous buffer of choice. VDM11 has a solubility of approximately 0.25 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Numerous analogs of arachidonoyl ethanolamide (AEA) potentiate its biological activity. This potentiation is ascribed either to inhibition of AEA reuptake into neurons, or inhibition of fatty amide acyl hydrolase (FAAH) within the neurons.<sup>1,2</sup> VDM11 is an AEA transport inhibitor with essentially no activity on either the central cannabinoid receptor (CB<sub>1</sub>), peripheral cannabinoid receptor (CB<sub>2</sub>), or the vanilloid receptor 1 (VR<sub>1</sub>).<sup>3</sup> However, VDM11 inhibits FAAH and monoacylglycerol lipase (MAGL) and may act as an alternative FAAH substrate.<sup>4</sup> At a concentration of 3 μM, VDM11, like AM404, inhibits glutamergic synaptic transmission between hippocampal neurons.<sup>5</sup> The mechanism of this effect may be a direct action on sodium channels. Thus, the use of anandamide analogs as uptake inhibitors and interpretation of the results must be undertaken with care.

### References

1. Khanolkar, A.D. and Makriyannis, A. Structure-activity relationships of anandamide, an endogenous cannabinoid ligand. *Life Sci.* **65**, 607-616 (1999).
2. Deutsch, D.G., Glaser, S.T., Howell, J.M., *et al.* The cellular uptake of anandamide is coupled to its breakdown by fatty-acid amide hydrolase. *J. Biol. Chem.* **276**(10), 6967-6973 (2001).
3. De Petrocellis, L., Bisogno, T., Davis, B.J., *et al.* Overlap between the ligand recognition properties of the anandamide transporter and the VR1 vanilloid receptor: Inhibitors of anandamide uptake with negligible capsaicin-like activity. *FEBS* **483**, 52-56 (2000).
4. Vandervoort, S. and Fowler, C.J. Inhibition of fatty acid amide hydrolase and monoacylglycerol lipase by the anandamide uptake inhibitor VDM11: Evidence that VDM11 acts as an FAAH substrate. *Br. J. Pharmacol.* **145**, 885-893 (2005).
5. Kelley, B.G. and Thayer, S.A. Anandamide transport inhibitor AM404 and structurally related compounds inhibit synaptic transmission between rat hippocampal neurons in culture independent of cannabinoid CB<sub>1</sub> receptors. *Eur. J. Pharmacol.* **496**, 33-39 (2004).

#### 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, 03/07/2024

#### 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