

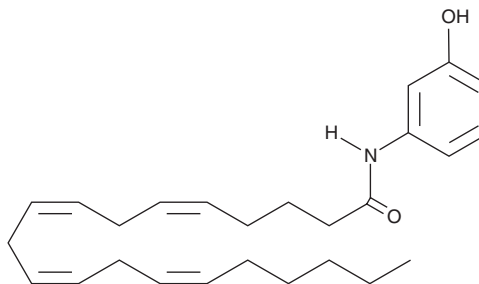
PRODUCT INFORMATION



N-(3-hydroxyphenyl)-Arachidonoyl amide

Item No. 10007704

CAS Registry No.: 183718-75-4
Formal Name: N-(3-hydroxyphenyl)-5Z,8Z,11Z,14Z-eicosatetraenamide
Synonyms: 3-HPA, 3-HPAA
MF: C₂₆H₃₇NO₂
FW: 395.6
Purity: ≥98%
UV/Vis.: λ_{max}: 246, 282 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

N-(3-hydroxyphenyl)-Arachidonoyl amide (3-HPA) 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 3-HPA in these solvents is approximately 20 mg/ml.

3-HPA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 3-HPA should be diluted with the aqueous buffer of choice. 3-HPA has a solubility of 0.3 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

3-HPA is an analog of AM404 (N-(4-hydroxyphenyl)-arachidonoyl amide), which is a selective inhibitor of carrier-mediated transport of arachidonoyl ethanolamide (AEA).¹ 3-HPA is metabolized by both cyclooxygenase-1 (COX-1) and COX-2 and was found to selectively and irreversibly inhibit COX-2 with an IC₅₀ value of 2 μM.²

References

1. Beltramo, M., Stella, N., Calignano, A., *et al.* Functional role of high-affinity anandamide transport, as revealed by selective inhibition. *Science* **277**(5329), 1094-1097 (1997).
2. Turman, M.V., Kinsgley, P.J., and Marnet, L.J. 9th International conference on eicosanoids and other bioactive lipids in cancer, inflammation, and related diseases. (2005).

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, 03/06/2024

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