

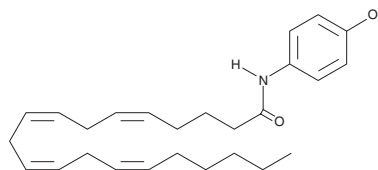
PRODUCT INFORMATION



AM404

Item No. 90060

CAS Registry No.: 183718-77-6
Formal Name: N-(4-hydroxyphenyl)-5Z,8Z,11Z,14Z-eicosatetrenamide
Synonyms: 4-HPA, N-(4-hydroxyphenyl)-Arachidonoyl Amide
MF: C₂₆H₃₇NO₂
FW: 395.6
Purity: ≥98%
UV/Vis.: λ_{max}: 250 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

AM404 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 AM404 in these solvents is approximately 50 mg/ml.

AM404 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of AM404 should be diluted with the aqueous buffer of choice. AM404 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AM404 is an analog of arachidonoyl ethanolamide (AEA) which potentiates the activity of endogenous AEA by blocking its re-uptake into presynaptic neurons. AM404 selectively inhibits the carrier-mediated transport of AEA without affecting anandamide hydrolysis. It inhibits the transport of AEA with an IC₅₀ of 1 μM in rat neurons and 5 μM in rat astrocytes. In *in vivo* models, AM404 enhances and prolongs exogenous AEA-induced analgesia at a dose of 10 mg/kg.¹

Reference

1. Beltramo, M., Stella, N., Calignano, A., *et al.* Functional role of high-affinity anandamide transport, as revealed by selective inhibition. *Science* **277**(5329), 1094-1096 (1997).

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, 02/28/2024

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