

Product Information



Arachidonoyl Glycine-d₈

Catalog No. 10007531

Formal Name: N-[1-oxo-5Z,8Z,11Z,14Z-eicosatetraenyl]-glycine,5,6,8,9,11,12,14,15-d₈

Synonyms: N-Arachidonoyl Glycine-d₈; NAGly-d₈

MF: C₂₂H₂₇D₈NO₃

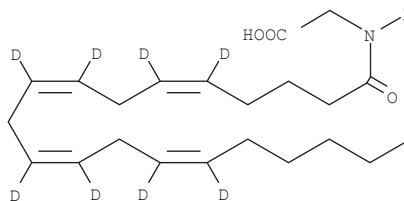
FW: 369.6

Chemical Purity: ≥97%

Deuterium Incorporation: ≤1% d₀

Stability: ≥1 year at -20°C

Supplied as: A solution in ethanol



Laboratory Procedures

Arachidonoyl glycine-d₈ (NAGly-d₈) contains eight deuterium atoms at the 5, 6, 8, 9, 11, 12, 14, and 15 positions. It is intended for use as an internal standard for the quantification of NAGly by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that NAGly-d₈ be stored as supplied at -20°C. It will be stable for at least two years.

NAGly-d₈ 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 NAGly-d₈ in these solvents is approximately 15 mg/ml.

NAGly-d₈ is used as an internal standard for the quantification of NAGly by stable isotope dilution MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

NAGly has been isolated from cell cultures treated with arachidonoyl ethanolamide (AEA),¹ from extracts of mammalian brain,^{2,3} and has also been synthesized as an analog of AEA for structure/activity testing.⁴ NAGly may be produced endogenously *via* oxidation of AEA or by transacylation of arachidonoyl coenzyme A. NAGly is reported to have analgesic activities in whole animal experiments.¹⁻³ Since it seems to be a very poor ligand for the CB₁ receptor,⁴ these effects are probably mediated *via* other signalling pathways.

References

1. Burstein, S.H., Rossetti, R.G., Yagen, B., *et al.* Oxidative metabolism of anandamide. *Prostaglandins and Other Lipid Mediators* **61**, 29-41 (2000).
2. Huang, S.M., Bisogno, T., Petros, T.J., *et al.* Identification and characterization of an endogenous anandamide-like compound: N-arachidonoylglycine (NAGly). *ICRS 2001 Symposium on the Cannabinoids* 78 (2001).
3. Huang, S.M., Bisogno, T., Petros, T.J., *et al.* Identification of a new class of molecules, the arachidonoyl amino acids, and characterization of one member that inhibits pain. *J. Biol. Chem.* **276**(46), 42639-42644 (2001).
4. Sheskin, T., Hanus, L., Slager, J., *et al.* Structural requirements for binding of anandamide-type compounds to the brain cannabinoid receptor. *J. Med. Chem.* **40**, 659-667 (1997).

Related Products

Arachidonoyl Ethanolamide - Cat. No. 90050 • Arachidonoyl Glycine - Cat. No. 90051 • N-Oleoylglycine - Cat. No. 90269 • Arachidonoyl Ethanolamide-d₈ - Cat. No. 390050

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our **website and in our catalog**.

Copyright Cayman Chemical Company, 08/21/2007