Product Information



N-Arachidonoyl Dopamine-d_o

1150000 (2 5

Item No. 10007431

CAC D . N

CAS Registry No.:	11)9908-42-3	
Formal Name:	N-[2-(3,4-dihydroxyphenyl)ethyl]-	
	5Z,8Z,11Z,14Z-eicosatetraenamide-	,OH
	5,6,8,9,11,12,14,15-d ₈	
Synonym:	NADA-d ₈	H, , , , ,
MF:	$C_{28}H_{33}D_8NO_3$	
FW:	447.6	
Chemical Purity:	≥98%	$\langle \lor \lor \lor \circ$
Deuterium		$\backslash - \land - \land \land \land$
Incorporation:	\geq 99% deuterated forms (d ₁ -d ₈); \leq 1% d ₀	
Stability:	≥1 year at -20°C	D D D
Supplied as:	A solution in ethanol	
UV/Vis.:	λ _{max} : 205, 283 nm	

Laboratory Procedures

N-Arachidonoyl dopamine-d₈ (NADA-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 NADA by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that NADA-d₈ be stored as supplied at -20°C. It should be stable for at least one year.

NADA-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. NADA-d₈ is miscible in these solvents.

NADA-d₈ is used as an internal standard for the quantification of NADA 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).

Several different arachidonoyl amino acids, including NADA, have been isolated and characterized from bovine brain.¹ NADA is the amide of the neurotransmitter dopamine and arachidonic acid. NADA is a CB1-selective cannabinoid agonist, inducing the typical tetrad of hypothermia, analgesia, catalepsy, and hypomotility in rats which exceeds that of anandamide (AEA).² NADA is a full agonist at the vanilloid receptor 1, but is inactive on the dopaminergic D1 and D2 receptors. NADA is also a potent inhibitor (IC₅₀ = $0.25 \ \mu$ M) of the proliferation of MCF-7 breast carcinoma cells. Recent reports of NADA's endothelium-dependent vasodilation indicate that some of its cannabinergic activities antagonized by SR141716A may be non-CB₁/CB₂ dependent.³

References

- 1. Huang, S.M., Bisogno, T., Petros, T.J., et al. Identification of a new class of molecules, the arachidonyl amino acids, and characterization of one member that inhibits pain. J. Biol. Chem. 276(46), 42639-42644 (2001).
- 2. Bezuglov, V., Bobrov, M., Gretskaya, N., et al. Synthesis and biological evaluation of novel amides of polyunsaturated fatty acids with dopamine. Bioorg. Medicinal Chem. Letters 11, 447-449 (2001).
- 3. Randall, M. and Maxey, K.M. Personal Communication.

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10007431

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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 Safety Data Sheet, which has been sent via email to 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

purpose, suitability and interchantability which extends beyond the description of the chemicals interior. Cayman warrants only to the original customer that the interchantability matches 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, indirectal 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. Buyers' **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a <u>refund</u> of the purchase price, or at Cayman's option, the <u>replacement</u>, 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 tis destination. Failure of Buyer to give said notice within this result of the purchase price are or a cit or travel or travel or the second price of the purchase price are and it in the second price of the purchase price are arrival of the material at its destination. Failure of Buyer to give said notice within this result of the purchase price are arrival of the material at its destination. Failure of Buyer to give said notice within this result of the purchase price are arrival of the material at its destination. Failure of Buyer to give said notice within this result of the purchase price are arrival of the material dat is destination. Failure of Buyer to give said notice within this result of the purchase price.

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, 11/19/2013

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