

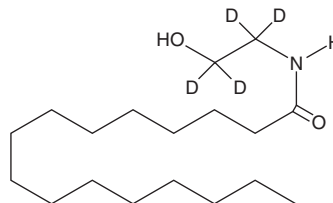
# Product Information



## Palmitoyl Ethanolamide-d<sub>4</sub>

Item No. 9000551

<b>Formal Name:</b>	N-(2-hydroxyethyl-1,1,2,2-d <sub>4</sub> )-hexadecanamide
<b>Synonyms:</b>	Palmidrol-d <sub>4</sub> , PEA-hydroxyethyl-1,1,2,2-d <sub>4</sub>
<b>MF:</b>	C <sub>18</sub> H <sub>33</sub> D <sub>4</sub> NO <sub>2</sub>
<b>FW:</b>	303.5
<b>Chemical Purity:</b>	≥98%
<b>Deuterium Incorporation:</b>	≤1% d <sub>0</sub>
<b>Stability:</b>	≥1 years at -20°C
<b>Supplied as:</b>	A solution in ethanol



### Laboratory Procedures

Palmitoyl ethanolamide-d<sub>4</sub> (PEA-d<sub>4</sub>) contains four deuterium atoms at the hydroxyethyl 1, 1', 2, and 2' positions. It is intended for use as an internal standard for the quantification of PEA by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that PEA-d<sub>4</sub> be stored as supplied at -20°C. It should be stable for at least one year.

PEA-d<sub>4</sub> 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 (DMF) purged with an inert gas can be used. The solubility of PEA-d<sub>4</sub> in DMSO is approximately 5 mg/ml and approximately 10 mg/ml in DMF.

PEA-d<sub>4</sub> is used as an internal standard for the quantification of PEA 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).

Palmitoyl ethanolamide (PEA) is an endogenous cannabinoid found in brain, liver, and other mammalian tissues.<sup>1</sup> PEA has also been isolated from egg yolk, and found to have anti-anaphylactic and anti-inflammatory activity *in vitro*.<sup>2</sup> PEA is an endocannabinoid which has been shown to significantly elevate cAMP in cells expressing CB<sub>2</sub> receptors. However, its affinity for CB<sub>2</sub> receptors is relatively low, at about 10 μM. CB<sub>1</sub> receptors have no appreciable affinity for PEA.<sup>3</sup>

### References

1. Bachur, N.R., Masek, K., Melmon, K.L., *et al.* Fatty acid amides of ethanolamine in mammalian tissues. *J. Biol. Chem.* **240**, 1019-1024 (1965).
2. Ganley, O.H., Graessle, O.E., Robinson, H.J., *et al.* Anti-inflammatory activity of compounds obtained from egg yolk, peanut oil, and soybean lecithin. *J. Lab. Clin. Med.* **51**, 709-714 (1958).
3. Devane, W.A., Hanus, L., Breuer, A., *et al.* Isolation and structure of a brain constituent that binds to the cannabinoid receptor. *Science* **258**, 1946-1949 (1992).

### Related Products

Prostaglandin D<sub>2</sub> Ethanolamide - Item No. 12012 • Prostaglandin E<sub>1</sub> Ethanolamide - Item No. 13012 • Prostaglandin E<sub>2</sub> Ethanolamide - Item No. 14012 • Arachidonyl Ethanolamide - Item No. 90050 • R-1 Methanandamide - Item No. 90070 • S-1 Methanandamide - Item No. 90072 • R-2 Methanandamide - Item No. 90074 • S-2 Methanandamide - Item No. 90076 • Linoleyl Ethanolamide - Item No. 90155 • Mead Acid Ethanolamide - Item No. 90195 • α-Linolenyl Ethanolamide - Item No. 90215 • Dihomo-γ-Linolenyl Ethanolamide - Item No. 90235 • Stearoyl Ethanolamide - Item No. 90245 • Oleoyl Ethanolamide - Item No. 90265 • Palmitoyl Ethanolamide - Item No. 90350 • Docosotetraenyl Ethanolamide - Item No. 90385 • Palmitoyl Ethanolamide-d<sub>4</sub> - Item No. 10007824

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

#### 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 *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 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, 08/24/2010

### 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