

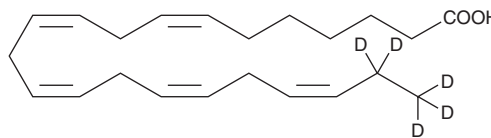
PRODUCT INFORMATION



Docosapentaenoic Acid-d₅

Item No. 20913

CAS Registry No.: 1215168-70-9
Formal Name: 7Z,10Z,13Z,16Z,19Z-docosapentaenoic-21,21,22,22,22-d₅ acid
Synonyms: DPA-d₅, n-3 DPA-d₅, FA 22:5-d₅
MF: C₂₂H₂₉D₅O₂
FW: 335.5
Chemical Purity: ≥98% (Docosapentaenoic Acid)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
Supplied as: A 500 µg/ml 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

Docosapentaenoic acid-d₅ (DPA-d₅) is intended for use as an internal standard for the quantification of DPA (Item No. 90165) by GC- or LC-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).

DPA-d₅ 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 DPA-d₅ in these solvents is approximately 100 mg/ml. The solubility of DPA-d₅ in 0.1 M Na₂CO₃ is approximately 1 mg/ml. DPA-d₅ is miscible in ethanol.

Description

DPA is an ω-3 fatty acid found in fish oils. It is a minor constituent of the total serum unsaturated fatty acids in humans, ranging from 0.1 to 1%, and increases on dietary supplementation.¹

Reference

1. Marckmann, P., Lassen, A., Haraldsdóttir, J., et al. Biomarkers of habitual fish intake in adipose tissue. *Am. J. Clin. Nutr.* **62**(5), 956-959 (1995).

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, 07/29/2025

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