

PRODUCT INFORMATION



Docosapentaenoic Acid (sodium salt)

Item No. 21907

Formal Name: 7Z,10Z,13Z,16Z,19Z-docosapentaenoic acid, monosodium salt

Synonyms: DPA, n-3 DPA, FA 22:5

MF: $C_{22}H_{33}O_2 \cdot Na$

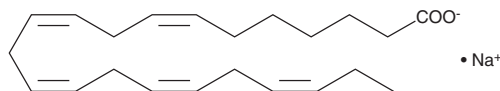
FW: 352.5

Purity: $\geq 98\%$

Supplied as: A crystalline solid

Storage: $-80^{\circ}C$

Stability: ≥ 1 year



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

Laboratory Procedures

Docosapentaenoic acid (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the docosapentaenoic acid (sodium salt) in the solvent of choice, which should be purged with an inert gas. Docosapentaenoic acid (sodium salt) is soluble in the organic solvent ethanol at a concentration of approximately 1.5 mg/ml.

Docosapentaenoic acid (sodium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, docosapentaenoic acid (sodium salt) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Docosapentaenoic acid (sodium salt) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Docosapentaenoic acid 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, 04/03/2024

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