

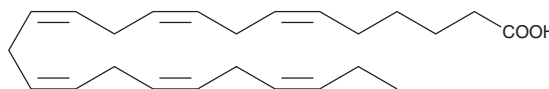
# PRODUCT INFORMATION



## 6(Z),9(Z),12(Z),15(Z),18(Z),21(Z)-Tetracosahexaenoic Acid

Item No. 10005165

**CAS Registry No.:** 68378-49-4  
**Formal Name:** 6Z,9Z,12Z,15Z,18Z,21Z-tetracosahexaenoic acid  
**Synonyms:** 24:6 (n-3), FA 24:6, Nisinic Acid, all-cis-6,9,12,15,18,21-Tetracosahexaenoic Acid  
**MF:** C<sub>24</sub>H<sub>36</sub>O<sub>2</sub>  
**FW:** 356.5  
**Purity:** ≥95%  
**Supplied as:** A solution in methyl acetate  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

6(Z),9(Z),12(Z),15(Z),18(Z),21(Z)-Tetracosahexaenoic acid is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the 6(Z),9(Z),12(Z),15(Z),18(Z),21(Z)-tetracosahexaenoic acid under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 6(Z),9(Z),12(Z),15(Z),18(Z),21(Z)-tetracosahexaenoic acid in these solvents is approximately 50 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of 6(Z),9(Z),12(Z),15(Z),18(Z),21(Z)-tetracosahexaenoic acid is needed, it can be prepared by evaporating the 6(Z),9(Z),12(Z),15(Z),18(Z),21(Z)-tetracosahexaenoic acid and directly dissolving the neat oil in aqueous buffers. The solubility of 6(Z),9(Z),12(Z),15(Z),18(Z),21(Z)-tetracosahexaenoic acid in PBS, pH 7.2, is approximately 100 µg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

6(Z),9(Z),12(Z),15(Z),18(Z),21(Z)-Tetracosahexaenoic acid is a very long chain polyunsaturated fatty acid (VLCPUFA) that has been found in tunny, herring, cod-liver, pilot whale, sardine, and squalus-sucii-liver oils, bovine retina, and as a component of triglycerides and cholesterol esters in mouse and rat testis.<sup>1-3</sup> It is produced from dietary linolenic acid via a series of elongation and Δ<sup>6</sup>-desaturation reactions in fish and rat brain and undergoes β-oxidation to form docosahexaenoic acid (DHA; Item No. 90310).<sup>4,5</sup>

### References

1. Ueno, S.-I. and Yonese, C. On the occurrence of new highly unsaturated fatty acids, C<sub>26</sub>H<sub>40</sub>O<sub>2</sub> and C<sub>26</sub>H<sub>42</sub>O<sub>2</sub>, in tunny oil. *J. Soc. Chem. Ind.* **11(7)**, 437-442 (1936).
2. Rötstein, N.P. and Aveldano, M.I. Synthesis of very long chain (up to 36 carbon) tetra, penta and hexaenoic fatty acids in retina. *Biochem. J.* **249(1)**, 191-200 (1988).
3. Furland, N.E., Maldonado, and Aveldano, M.I. Very long chain PUFA in murine testicular triglycerides and cholesterol esters. *Lipids* **38(1)**, 73-80 (2003).
4. Williard, D.E., Harmon, S.D., Kaduce, T.L., et al. Docosahexaenoic acid synthesis from n-3 polyunsaturated fatty acids in differentiated rat brain astrocytes. *J. Lipid Res.* **42(9)**, 1368-76 (2001).
5. Mourente, G. and Tocher, D.R. *In vivo* metabolism of [1-<sup>14</sup>C]linolenic acid (18:3(n-3)) and [1-<sup>14</sup>C]eicosapentaenoic acid (20:5(n-3)) in a marine fish: Time-course of the desaturation/elongation pathway. *Biochim. Biophys. Acta.* **1212(1)**, 109-118 (1994).

#### 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, 03/19/2024

#### 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