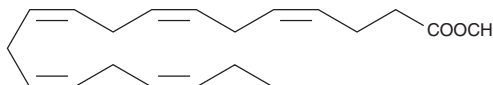


PRODUCT INFORMATION

4(Z),7(Z),10(Z),13(Z),16(Z)-Nonadecapentaenoic Acid methyl ester

Item No. 19472

CAS Registry No.: 2055816-84-5
Formal Name: 4(Z),7(Z),10(Z),13(Z),16(Z)-nonadecapentaenoic acid, methyl ester
Synonyms: C20:5 (cis-4,7,10,13,16) methyl ester, SFE 20:5
MF: C₂₀H₃₀O₂
FW: 302.5
Purity: ≥95%
Supplied as: A 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

4(Z),7(Z),10(Z),13(Z),16(Z)-Nonadecapentaenoic acid methyl ester 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 ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 4(Z),7(Z),10(Z),13(Z),16(Z)-nonadecapentaenoic acid methyl ester in ethanol is approximately 500 mg/ml and approximately 100 mg/ml in DMSO and DMF.

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 4(Z),7(Z),10(Z),13(Z),16(Z)-nonadecapentaenoic acid methyl ester is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 4(Z),7(Z),10(Z),13(Z),16(Z)-nonadecapentaenoic acid methyl ester in PBS, pH 7.2, is approximately 0.15 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

4(Z),7(Z),10(Z),13(Z),16(Z)-Nonadecapentaenoic acid methyl ester is an esterified form of 4(Z),7(Z),10(Z),13(Z),16(Z)-nonadecapentaenoic acid (Item No. 19471), which is a byproduct in the synthesis of 4(Z),7(Z),10(Z),13(Z)-hexadecatetraenoic acid methyl ester (Item No. 9002661).

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, 02/14/2024

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