

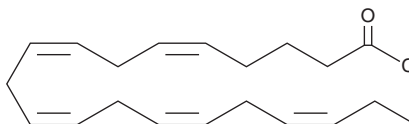
PRODUCT INFORMATION



Eicosapentaenoyl Chloride

Item No. 22721

CAS Registry No.: 98770-65-1
Formal Name: (5Z,8Z,11Z,14Z,17Z)-5,8,11,14,17-eicosapentaenoyl chloride
Synonyms: Eicosapentaenoic Acid Chloride, EPA Chloride
MF: C₂₀H₂₉ClO
FW: 320.9
Purity: ≥95%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Eicosapentaenoyl chloride is supplied as a liquid. A stock solution may be made by dissolving the eicosapentaenoyl chloride in the solvent of choice, which should be purged with an inert gas. Eicosapentaenoyl chloride is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of eicosapentaenoyl chloride in these solvents is approximately 10 mg/ml.

Description

Eicosapentaenoyl chloride is a derivative of eicosapentaenoic acid (Item Nos. 90110 | 90110.1 | 21908). It has been used in the synthesis of fatty acid conjugates to enhance lipophilicity and cell permeability of bioactive compounds such as (-)-epigallocatechin gallate (EGCG; Item No. 70935) and salicylic acid.^{1,2}

References

1. Zhong, Y., Chiou, Y.-S., Pan, M.-H., *et al.* Anti-inflammatory activity of lipophilic epigallocatechin gallate (EGCG) derivatives in LPS-stimulated murine macrophages. *Food Chem.* **134**(2), 742-748 (2012).
2. Vu, C.-B., Bemis, J.E., Benson, E., *et al.* Synthesis and characterization of fatty acid conjugates of niacin and salicylic acid. *J. Med. Chem.* **59**(3), 1217-1231 (2016).

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, 04/06/2020

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