

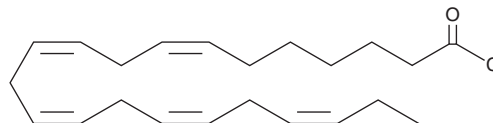
# PRODUCT INFORMATION



## Docosapentaenoyl Chloride

Item No. 22720

CAS Registry No.: 960593-70-8  
Formal Name: 7Z,10Z,13Z,16Z,19Z-  
docosapentaenoyl chloride  
MF:  $C_{22}H_{33}ClO$   
FW: 349.0  
Purity:  $\geq 95\%$   
Supplied as: A neat oil  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



**Special Conditions:** Do not store in ethanol, product will esterify.

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

### Laboratory Procedures

Docosapentaenoyl chloride is supplied as a neat oil. A stock solution may be made by dissolving the docosapentaenoyl chloride in the solvent of choice, which should be purged with an inert gas. Docosapentaenoyl chloride is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of docosapentaenoyl chloride in these solvents is approximately 15 and 20 mg/ml, respectively.

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. Organic solvent-free aqueous solutions of docosapentaenoyl chloride can be prepared by directly dissolving the neat oil in aqueous buffers. The solubility of docosapentaenoyl chloride in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Docosapentaenoyl chloride is a derivative of the fatty acid docosapentaenoic acid (DPA; Item No. 90165 | 21907), an  $\omega$ -3 fatty acid found in fish oils, that is more reactive than the free fatty acid.

#### 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, 12/16/2022

#### 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