

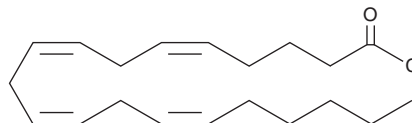
# PRODUCT INFORMATION



## Arachidonoyl Chloride

Item No. 18078

**CAS Registry No.:** 57303-04-5  
**Formal Name:** 5Z,8Z,11Z,14Z-eicosatetraenoyl chloride  
**Synonym:** Arachidonic Acid Chloride  
**MF:** C<sub>20</sub>H<sub>31</sub>ClO  
**FW:** 322.9  
**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

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

### Description

Arachidonoyl chloride is a derivative of arachidonic acid (Item No. 90010). It has been used as an intermediate in the synthesis of arachidonic acid derivatives.<sup>1</sup>

### Reference

1. Clapper, J.R., Duranti, A., Tontini, A., *et al.* The fatty-acid amide hydrolase inhibitor URB597 does not affect triacylglycerol hydrolysis in rat tissues. *Pharmacol. Res.* **54(5)**, 341-344 (2006).

#### 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, 10/10/2022

#### 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