

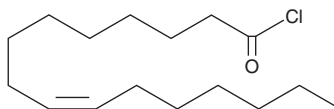
# PRODUCT INFORMATION



## Palmitoleoyl Chloride

Item No. 22723

CAS Registry No.: 40426-22-0  
Formal Name: 9Z-hexadecenoyl chloride  
Synonym: Palmitoleic Acid Chloride  
MF:  $C_{16}H_{29}ClO$   
FW: 272.9  
Purity:  $\geq 98\%$   
Supplied as: A liquid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 2$  years



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

### Laboratory Procedures

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

### Description

Palmitoleoyl chloride is a derivative of palmitoleic acid (Item Nos. 10009871 | 21911). It has been used as an intermediate in the synthesis of a variety of compounds, including antiproliferative and anti-inflammatory compounds.<sup>1,2</sup>

### References

1. Hassan, A.H.E., Park, H.R., Yoon, Y.M., *et al.* Antiproliferative 3-deoxysphingomyelin analogs: Design, synthesis, biological evaluation and molecular docking of pyrrolidine-based 3-deoxysphingomyelin analogs as anticancer agents. *Bioorg. Chem.* **84**, 444-455 (2019).
2. Fox, D.J., Reckless, J., Wilbert, S.M., *et al.* Identification of 3-(acylamino)azepan-2-ones as stable broad-spectrum chemokine inhibitors resistant to metabolism in vivo. *J. Med. Chem.* **48**(3), 867-874 (2005).

#### 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, 04/19/2019

#### 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