

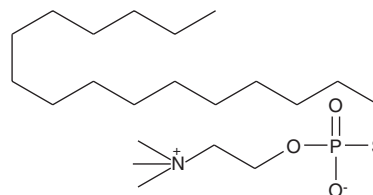
# PRODUCT INFORMATION



## thio-Miltefosine

Item No. 10009813

CAS Registry No.: 943022-11-5  
Formal Name: 1-hexadecyl-thio-phosphorylcholine  
MF:  $C_{21}H_{46}NO_3PS$   
FW: 423.6  
Purity:  $\geq 98\%$   
Supplied as: A neat oil  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 1$  year



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

### Laboratory Procedures

thio-Miltefosine is supplied as a neat oil. A stock solution may be made by dissolving the thio-miltefosine in the solvent of choice, which should be purged with an inert gas. thio-Miltefosine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of thio-miltefosine in these solvents is approximately 1, 0.5, and 0.05 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 thio-miltefosine can be prepared by directly dissolving the neat oil in aqueous buffers. The solubility of thio-miltefosine in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

thio-Miltefosine is a sulfur-containing derivative of miltefosine (Item No. 63280).<sup>1</sup>

### Reference

1. Perrissoud, D., Pietras, M., and Engel, J. Novel alkyl phospholipid derivatives with reduced cytotoxicity and uses thereof. *Zentaris GmbH. WO 2007/071658 A2* (2007).

#### 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, 06/30/2021

#### 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