

PRODUCT INFORMATION



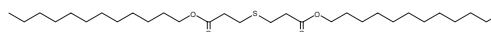
Dilauryl thiodipropionate

Item No. 42578

CAS Registry No.: 123-28-4

Formal Name: 3,3'-thiobis-propanoic acid,
1,1'-didodecyl ester

Synonyms: Didodecyl 3,3'-Thiodipropionic Acid,
Dilauryl 3,3'-Thiodipropanoic Acid,
Dilauryl β,β' -Thiodipropionic Acid,
DLTDP, DLTP, NSC 65494



MF: $C_{30}H_{58}O_4S$

FW: 514.8

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

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

Laboratory Procedures

Dilauryl thiodipropionate is supplied as a solid. A stock solution may be made by dissolving the dilauryl thiodipropionate in the solvent of choice, which should be purged with an inert gas. Dilauryl thiodipropionate is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

Dilauryl thiodipropionate is an antioxidant.¹ It inhibits performic acid-induced oxidation of ethyl oleate, nonanal, and cholesterol. Formulations containing dilauryl thiodipropionate have been used as antioxidants in cosmetics.

Reference

1. Karahadian, C. and Lindsay, R.C. Evaluation of the mechanism of dilauryl thiodipropionate antioxidant activity. *J. Am. Oil Chem. Soc.* **65**, 1159-1165 (2014).

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, 01/29/2025

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