

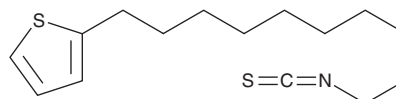
PRODUCT INFORMATION



Thienyldecyl isothiocyanate

Item No. 19653

CAS Registry No.: 288323-41-1
Formal Name: 2-(10-isothiocyanatodecyl)-thiophene
MF: $C_{15}H_{23}NS_2$
FW: 281.5
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 235 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

Thienyldecyl isothiocyanate is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of thienyldecyl isothiocyanate in ethanol is approximately 30 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Thienyldecyl isothiocyanate is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Thienyldecyl isothiocyanate is an analog of thienylbutyl isothiocyanate that, like many isothiocyanates, demonstrates antiproliferative activity against cancer cells presumably by modulation of xenobiotic-metabolizing enzymes, such as by inhibition of cytochrome P450, and/or by induction of phase II detoxifying enzymes.¹

Reference

1. Yu, R., Mandlekar, S., Harvey, K.J., *et al.* Chemopreventive isothiocyanates induce apoptosis and caspase-3-like protease activity. *Cancer Res.* **58(3)**, 402-408 (1998).

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, 12/15/2022

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