

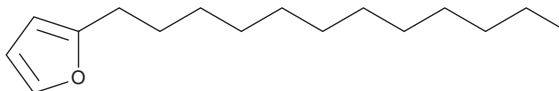
PRODUCT INFORMATION



2-n-Dodecylfuran

Item No. 39269

CAS Registry No.: 75308-12-2
Formal Name: 2-dodecyl-furan
MF: $C_{16}H_{28}O$
FW: 236.4
Purity: $\geq 98\%$
UV/Vis.: $\lambda_{\max}$: 216 nm
Supplied as: A neat oil
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

2-n-Dodecylfuran is supplied as a neat oil. A stock solution may be made by dissolving the 2-n-dodecylfuran in the solvent of choice, which should be purged with an inert gas. 2-n-Dodecylfuran is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2-n-dodecylfuran in these solvents is approximately 25, 5, and 50 mg/ml, respectively.

Description

2-n-Dodecylfuran is an intermediate in the synthesis of thermally cleavable surfactants.¹

Reference

1. McElhanon, J.R., Zifer, T., Kline, S.R., *et al.* Thermally cleavable surfactants based on furan-maleimide Diels-Alder adducts. *Langmuir* **21**(8), 3259-3266 (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, 10/03/2023

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