

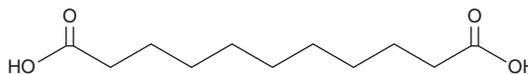
PRODUCT INFORMATION



Undecanedioic Acid

Item No. 36465

CAS Registry No.: 1852-04-6
Synonyms: FA 11:1;O2,
Hendecanedioic Acid,
NSC 400241, UDA,
1,9-Nonanedicarboxylic Acid,
1,11-Undecanedioic Acid
MF: C₁₁H₂₀O₄
FW: 216.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Undecanedioic acid is supplied as a solid. A stock solution may be made by dissolving the undecanedioic acid in the solvent of choice, which should be purged with an inert gas. Undecanedioic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of undecanedioic acid in DMF is approximately 16 mg/ml and approximately 20 mg/ml in ethanol and DMSO.

Description

Undecanedioic acid is a medium-chain dicarboxylic acid that has been found in *S. officinalis*.^{1,2} It is active against *T. rubrum*, *T. mentagrophytes*, and *M. canis* when used at a concentration of 1 mM.² Undecanedioic acid has been used in the synthesis of proteolysis-targeting chimeras (PROTACs) with anticancer activity in mutant EGFR-expressing cancer cells.³ It has been found in urban and industrial aerosols of fine particulate matter less than 2.5 μm (PM_{2.5}).⁴

References

1. Aldal'In, H.K., Wedian, F., Al-Mazaideh, G.M., *et al.* Comparative analysis of phytochemical composition of ethanolic extract of Jordanian *Silvia officinalis*. *Pak. J. Biol. Sci.* **23**(8), 989-994 (2022).
2. Brasch, J. and Friege, B. Dicarboxylic acids affect the growth of dermatophytes *in vitro*. *Acta Derm. Venereol.* **74**(5), 347-350 (1994).
3. Li, Q., Guo, Q., Wang, S., *et al.* Design and synthesis of proteolysis targeting chimeras (PROTACs) as an EGFR degrader based on CO-1686. *Eur. J. Med. Chem.* **238**, 114455 (2022).
4. Crenn, V., Fronval, I., Petitprez, D., *et al.* Fine particles sampled at an urban background site and an industrialized coastal site in Northern France - Part 1: Seasonal variations and chemical characterization. *Sci. Total Environ.* **578**, 203-218 (2017).

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, 02/20/2024

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