

PRODUCT INFORMATION



DTUN

Item No. 32742

Formal Name: (E)-1,2-bis((2-methyldecan-2-yl)oxy)diazene

MF: C₂₂H₄₆N₂O₂

FW: 370.6

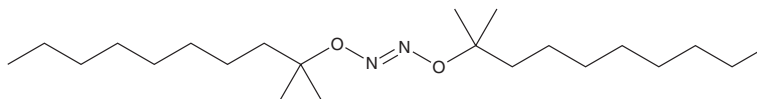
Purity: ≥95%

UV/Vis.: λ_{max}: 227 nm

Supplied as: A film

Storage: -80°C

Stability: ≥2 years



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

Laboratory Procedures

DTUN is supplied as a film. A stock solution may be made by dissolving the DTUN in the solvent of choice, which should be purged with an inert gas. DTUN is soluble in the organic solvent ethanol at a concentration of approximately 50 mg/ml.

Description

DTUN is a lipophilic hyponitrite radical initiator.¹ It initiates co-oxidation of STY-BODIPY-embedded egg phosphatidylcholine liposomes when used at a concentration of 0.2 mM. DTUN has been used in fluorescence-enabled inhibited autoxidation (FENIX) assays for the development of ferroptosis inhibitors and the study of antioxidants in lipid membranes.^{1,2}

References

1. Shah, R., Farmer, L.A., Zilka, O., *et al.* Beyond DPPH: Use of fluorescence-enabled inhibited autoxidation to predict oxidative cell death rescue. *Cell Chem. Biol.* **26(11)**, 1594-1607 (2019).
2. Soula, M., Weber, R.A., Zilka, O., *et al.* Metabolic determinants of cancer cell sensitivity to canonical ferroptosis inducers. *Nat. Chem. Biol.* **16(12)**, 1351-1360 (2020).

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, 06/23/2022

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