

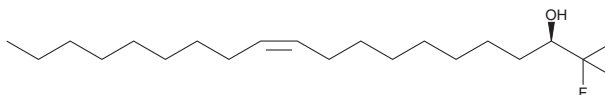
PRODUCT INFORMATION



1(R)-(Trifluoromethyl)oleyl alcohol

Item No. 31796

CAS Registry No.: 2644636-20-2
Formal Name: (2R,10Z)-1,1,1-trifluoro-10-nonadecen-2-ol
MF: C₁₉H₃₅F₃O
FW: 336.5
Purity: ≥95%
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.

Description

1(R)-(Trifluoromethyl)oleyl alcohol is an analog of oleic acid (Item Nos. 90260 | 24659).¹ It inhibits ferroptosis induced by erastin (Item No. 17754) in primary fibroblasts isolated from patients with Friedreich ataxia, a neuro- and cardiodegenerative disorder characterized by loss or impaired activity of frataxin (FXN), when used at concentrations of 5, 10, or 20 μM. 1(R)-(Trifluoromethyl)oleyl alcohol (5 μM) reduces lipid peroxidation induced by FXN siRNA knockdown in NBT human myoblasts.

Reference

1. Cotticelli, M.G., Forestieri, R., Xia, S., *et al.* Identification of a novel oleic acid analog with protective effects in multiple cellular models of Friedreich ataxia. *ACS Chem. Neurosci.* **11(17)**, 2535-2542 (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, 01/25/2024

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