

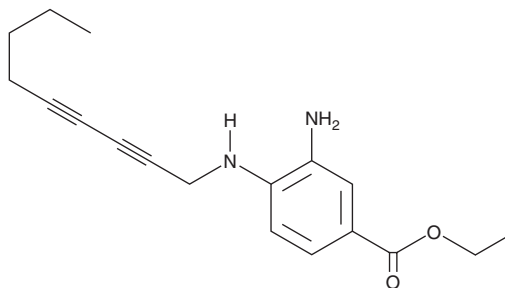
PRODUCT INFORMATION



Ferrostatin-1 Diyne

Item No. 25492

CAS Registry No.: 2226204-90-4
Formal Name: 3-amino-4-(2,4-nonadiyn-1-ylamino)-benzoic acid, ethyl ester
MF: C₁₈H₂₂N₂O₂
FW: 298.4
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 241, 290, 319 nm
Supplied as: A crystalline 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

Ferrostatin-1 diyne is supplied as a crystalline solid. A stock solution may be made by dissolving the ferrostatin-1 diyne in the solvent of choice, which should be purged with an inert gas. Ferrostatin-1 diyne is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ferrostatin-1 diyne in ethanol and DMSO is approximately 10 mg/ml and approximately 30 mg/ml in DMF.

Ferrostatin-1 diyne is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ferrostatin-1 diyne should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Ferrostatin-1 diyne has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ferrostatin-1 diyne is a ferroptosis inhibitor and analog of ferrostatin-1 (Item No. 17729).¹ It inhibits cell death induced by erastin (Item No. 17754) in HT-1080 cells when used at concentrations ranging from 0.01 to 1 μM. Ferrostatin-1 diyne contains a diyne small vibrational tag and has been used to assess localization of ferrostatin-1 in HT-1080 cells using stimulated Raman scattering microscopy.

Reference

1. Gaschler, M.M., Hu, F., Feng, H., *et al.* Determination of the subcellular localization and mechanism of action of ferrostatins in suppressing ferroptosis. *ACS Chem. Biol.* **13**(4), 1013-1020 (2018).

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, 02/16/2024

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