

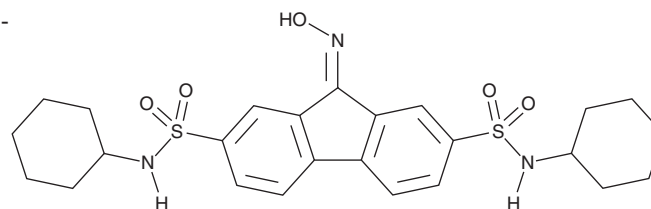
PRODUCT INFORMATION



FIN56

Item No. 25180

CAS Registry No.: 1083162-61-1
Formal Name: N²,N⁷-dicyclohexyl-9-(hydroxyimino)-9H-fluorene-2,7-disulfonamide
MF: C₂₅H₃₁N₃O₅S₂
FW: 517.7
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 267, 305 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

FIN56 is supplied as a crystalline solid. A stock solution may be made by dissolving the FIN56 in the solvent of choice, which should be purged with an inert gas. FIN56 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of FIN56 in these solvents is approximately 30 mg/ml.

Description

FIN56 is an inducer of ferroptosis and a derivative of CIL56 (Item No. 19287).¹ It selectively induces ferroptosis in BJ cells overexpressing oncogenic HRAS^{G12V} via acetyl-CoA carboxylase-dependent reduction in expression of GPX4 protein. This effect is reversed by the ferroptosis inhibitors deferoxamine (Item No. 14595) and α-tocopherol as well as overexpression of the fusion protein GFP-GPX4. FIN56 also binds to and activates squalene synthase in a GPX4 degradation-independent manner.

Reference

1. Shimada, K., Souta, R., Kaplan, A., *et al.* Global survey of cell death mechanisms reveals metabolic regulation of ferroptosis. *Nat. Chem. Biol.* **12**(7), 497-503 (2016).

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, 11/29/2022

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