

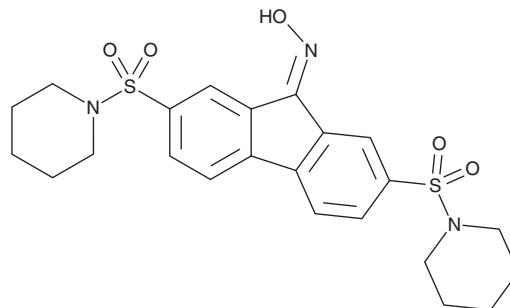
PRODUCT INFORMATION



CIL56

Item No. 19287

CAS Registry No.: 300802-28-2
Formal Name: 2,7-bis(1-piperidinylsulfonyl)-9H-fluoren-9-one, oxime
MF: C₂₃H₂₇N₃O₅S₂
FW: 489.6
Purity: ≥95%
UV/Vis.: λ_{max}: 224, 267, 308 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

CIL56 is supplied as a crystalline solid. A stock solution may be made by dissolving the CIL56 in the solvent of choice, which should be purged with an inert gas. CIL56 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CIL56 in ethanol is approximately 0.15 mg/ml and 10 mg/ml in DMSO and DMF.

CIL56 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CIL56 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CIL56 has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CIL56 is a small molecule that induces non-apoptotic cell death in a manner dependent on acetyl-CoA carboxylase 1 (ACC1), the rate-limiting enzyme in fatty acid synthesis.¹ HT-1080 cells lacking ACC1 exhibit 5-fold resistance to CIL56 treatment, indicating ACC1 activity sensitizes cells to CIL56-induced cell death. CIL56 (6.5 μM) induces accumulation of long-chain saturated, monounsaturated, and polyunsaturated fatty acids in HT-1080 cells *in vitro*.

Reference

1. Dixon, S.J., Winter, G.E., Musavi, L.S., *et al.* Human haploid cell genetics reveals roles for lipid metabolism genes in nonapoptotic cell death. *ACS Chem. Biol.* **10**(7), 1604-1609 (2015).

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