

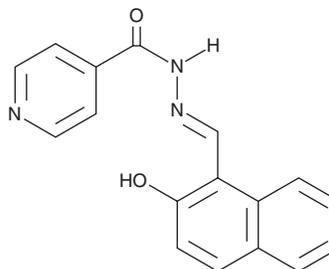
PRODUCT INFORMATION



AS-8351

Item No. 20811

CAS Registry No.: 796-42-9
Formal Name: 2-[(2-hydroxy-1-naphthalenyl)methylene]hydrazide, 4-pyridinecarboxylic acid
Synonyms: NCI 51355, NSC 51355
MF: C₁₇H₁₃N₃O₂
FW: 291.3
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 257, 327, 367 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

AS-8351 is supplied as a crystalline solid. A stock solution may be made by dissolving the AS-8351 in the solvent of choice, which should be purged with an inert gas. AS-8351 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AS-8351 in these solvents is approximately 20 and 25 mg/ml, respectively.

Description

AS-8351 is a histone demethylase inhibitor. It has been used in combination with CHIR99021 (Item No. 13122), A 83-01 (Item No. 9001799), BIX01294 (Item No. 13124), SC-1 (Item No. 10009557), Y-27632 (Item No. 10005583), OAC2 (Item No. 14103), SU 16f (Item No. 19555), and JNJ-10198409 (Item No. 10008131) to induce reprogramming of human fetal lung fibroblasts into functional cardiomyocytes.¹

Reference

1. Cao, N., Huang, Y., Zheng, J., *et al.* Conversion of human fibroblasts into functional cardiomyocytes by small molecules. *Science* **352**(6290), 1216-1220 (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, 12/19/2022

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