

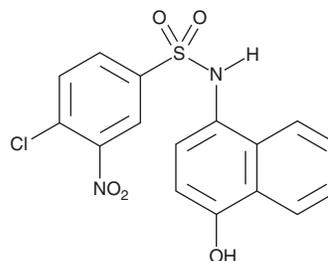
PRODUCT INFORMATION



SW155246

Item No. 21865

CAS Registry No.: 420092-79-1
Formal Name: 4-chloro-N-(4-hydroxy-1-naphthalenyl)-3-nitrobenzenesulfonamide
MF: C₁₆H₁₁ClN₂O₅S
FW: 378.8
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 304 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

SW155246 is supplied as a crystalline solid. A stock solution may be made by dissolving the SW155246 in the solvent of choice. SW155246 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of SW155246 in these solvents is approximately 30 mg/ml.

SW155246 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SW155246 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SW155246 has a solubility of approximately 0.16 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SW155246 is an inhibitor of DNA methyltransferase 1 (DNMT1; IC₅₀ = 1.2 μM).¹ It is selective for human DNMT1 over murine DNMT3 (IC₅₀ = 38 μM). SW155246 inhibits HeLa and A549 tumor cell growth in a redox-independent manner (IC₅₀s = 8-20 μM). It also inhibits genomic methylation in HeLa cells when used at concentrations ranging from 10 to 18.6 μM.

Reference

1. Kilgore, J.A., Du, X., Melito, L., *et al.* Identification of DNMT1 selective antagonists using a novel scintillation proximity assay. *J. Biol. Chem.* **288**(27), 19673-19684 (2013).

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, 11/14/2022

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