

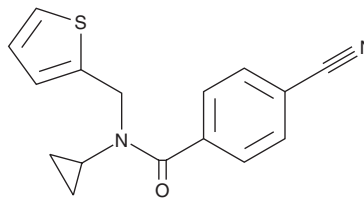
PRODUCT INFORMATION



MR837

Item No. 43711

CAS Registry No.: 1210906-48-1
Formal Name: 4-cyano-N-cyclopropyl-N-(2-thienylmethyl)-benzamide
MF: C₁₆H₁₄N₂OS
FW: 282.4
Purity: ≥98%
Supplied as: A 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

MR837 is supplied as a solid. A stock solution may be made by dissolving the MR837 in the solvent of choice, which should be purged with an inert gas. MR837 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

MR837 is an inhibitor of the histone-lysine N-methyltransferase NSD2 (IC₅₀ = 24.67 μM).¹ It binds to the NSD2 PWWP1 domain (K_d = 7 μM), which binds methylated histones, and inhibits the protein-protein interaction between NSD2 and histone H3.3 in U2OS cells (IC₅₀ = 17.3 μM).²

Reference

1. Li, N., Yang, H., Liu, K., *et al.* Structure-based discovery of a series of NSD2-PWWP1 inhibitors. *J. Med. Chem.* **65**(13), 9459-9477 (2022).
2. Ferreira de Freitas, R., Liu, Y., Szewczyk, M.M., *et al.* Discovery of small-molecule antagonists of the PWWP domain of NSD2. *J. Med. Chem.* **64**(3), 1584-1592 (2021).

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, 08/01/2025

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