

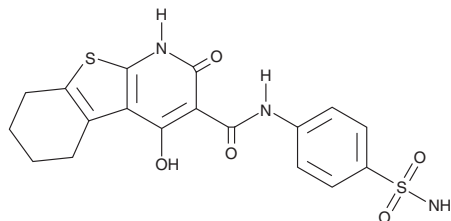
PRODUCT INFORMATION



M435-1279

Item No. 45401

CAS Registry No.: 1359431-16-5
Formal Name: N-[4-(aminosulfonyl)phenyl]-1,2,5,6,7,8-hexahydro-4-hydroxy-2-oxo-[1]benzothieno[2,3-b]pyridine-3-carboxamide
MF: C₁₈H₁₇N₃O₅S₂
FW: 419.5
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

M435-1279 is supplied as a solid. A stock solution may be made by dissolving the M435-1279 in the solvent of choice, which should be purged with an inert gas. M435-1279 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

M435-1279 is an inhibitor of ubiquitin-conjugating enzyme E2T (UBE2T).¹ It binds to UBE2T ($K_d = 50.5 \mu\text{M}$), inhibits UBE2T-mediated degradation of receptor for activated C kinase 1 (RACK1), and reduces hyperactive Wnt/ β -catenin signaling. M435-1279 reduces the viability of MKN45 and AGS gastric cancer cells when used at concentrations ranging from 4 to 32 μM . Intratumoral administration of M435-1279 (5 mg/kg per day) reduces tumor growth in an MKN45 mouse xenograft model.

Reference

1. Yu, Z., Jiang, X., Qin, L., *et al.* A novel UBE2T inhibitor suppresses Wnt/ β -catenin signaling hyperactivation and gastric cancer progression by blocking RACK1 ubiquitination. *Oncogene* **40(5)**, 1027-1042 (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, 04/15/2026

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