

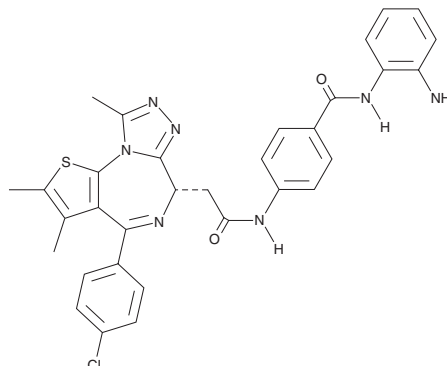
PRODUCT INFORMATION



TW9

Item No. 31588

CAS Registry No.: 2614417-90-0
Formal Name: (S)-N-(2-aminophenyl)-4-(2-(4-(4-chlorophenyl)-2,3,9-trimethyl-6H-thieno[3,2-f][1,2,4]triazolo[4,3-a][1,4]diazepin-6-yl)acetamido)benzamide
MF: C₃₂H₂₈ClN₇O₂S
FW: 610.1
Purity: ≥98%
UV/Vis.: λ_{max}: 259 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

TW9 is supplied as a crystalline solid. A stock solution may be made by dissolving the TW9 in the solvent of choice, which should be purged with an inert gas. TW9 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of TW9 in these solvents is approximately 10 and 20 mg/ml, respectively.

Description

TW9 is a dual inhibitor of bromodomain 2 (BD2) in bromodomain-containing protein 4 (BRD4) and histone deacetylase 1 (HDAC1; IC₅₀s = 0.074 and 0.29 μM, respectively).¹ It is selective for BD2 over BD1 in BRD4 (IC₅₀ = 0.72 μM) and for HDAC1 over HDAC2 (IC₅₀ = 2.5 μM). TW9 (50 nM) induces apoptosis in, and inhibits proliferation of, MIA PaCa-2 pancreatic cancer cells. It induces cell cycle arrest at the G₁ phase in HPAC pancreatic cancer cells when used at a concentration of 2 μM. TW9 acts synergistically with gemcitabine (Item No. 9003096) to reduce the viability of HPAC cells.

Reference

1. Zhang, X., Zegar, T., Weiser, T., *et al.* Characterization of a dual BET/HDAC inhibitor for treatment of pancreatic ductal adenocarcinoma. *Int. J. Cancer* **147**(10), 2847-2861 (2020).

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, 10/08/2021

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