

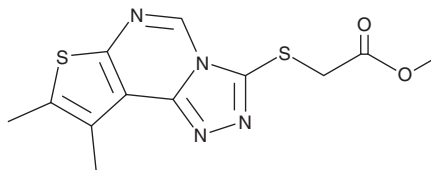
PRODUCT INFORMATION



BDW568

Item No. 39708

CAS Registry No.: 335401-44-0
Formal Name: 2-[(8,9-dimethylthieno[3,2-e]-1,2,4-triazolo[4,3-c]pyrimidin-3-yl)thio]-acetic acid, methyl ester
MF: C₁₂H₁₂N₄O₂S₂
FW: 308.4
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

BDW568 is supplied as a solid. A stock solution may be made by dissolving the BDW568 in the solvent of choice, which should be purged with an inert gas. BDW568 is soluble in chloroform, methanol, and acetonitrile.

Description

BDW568 is an agonist of stimulator of interferon genes (STING) and a prodrug form of BDW-OH.¹ It induces STING transcriptional activity in a reporter assay using THP-1 cells (EC₅₀ = 7.6 μM). BDW568 (50 μM) induces phosphorylation of TANK-binding kinase 1 (TBK1) and IFN regulatory factor 3 (IRF3) in THP-1 cells.

Reference

1. Zhao, J., Tang, Z., Selvaraju, M., et al. Cellular target deconvolution of small molecules using a selection-based genetic screening platform. *ACS Cent. Sci.* **8**(10), 1424-1434 (2022).

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, 11/26/2025

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