

PRODUCT INFORMATION

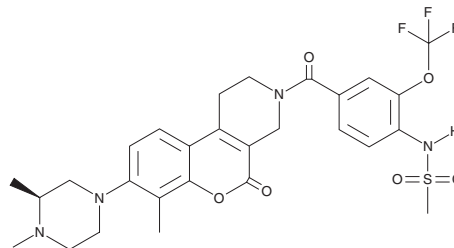


DS18561882

Item No. 33860

CAS Registry No.: 2227149-22-4
Formal Name: N-[4-[[8-[(3S)-3,4-dimethyl-1-piperazinyl]-1,5-dihydro-7-methyl-5-oxo-2H-[1]benzopyrano[3,4-c]pyridin-3(4H)-yl]carbonyl]-2-(trifluoromethoxy)phenyl]-methanesulfonamide

MF: C₂₈H₃₁F₃N₄O₆S
FW: 608.6
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

DS18561882 is supplied as a solid. A stock solution may be made by dissolving the DS18561882 in the solvent of choice, which should be purged with an inert gas. DS18561882 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

DS18561882 is an inhibitor of methylenetetrahydrofolate dehydrogenase 2 (MTHFD2; IC₅₀ = 0.0063 μM).¹ It is selective for MTHFD2 over MTHFD1 (IC₅₀ = 0.57 μM). DS18561882 inhibits the growth of MDA-MB-231 cells (GI₅₀ = 0.14 μM). *In vivo*, DS18561882 (100 and 300 mg/kg) reduces tumor growth in an MDA-MB-231 mouse xenograft model.

Reference

1. Kawai, J., Toki, T., Ota, M., *et al.* Discovery of a potent, selective, and orally available MTHFD2 inhibitor (DS18561882) with *in vivo* antitumor activity. *J. Med. Chem.* **62(22)**, 10204-10220 (2019).

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, 01/20/2025

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