

PRODUCT INFORMATION



DS44960156

Item No. 28366

CAS Registry No.: 2361327-08-2
Formal Name: 4-[(1,5-dihydro-5-oxo-2H-[1]benzopyrano[3,4-c]pyridin-3(4H)-yl)carbonyl]-benzoic acid

MF: C₂₀H₁₅NO₅

FW: 349.3

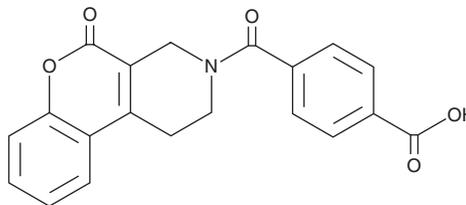
Purity: ≥98%

UV/Vis.: λ_{max}: 272 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

DS44960156 is supplied as a crystalline solid. A stock solution may be made by dissolving the DS44960156 in the solvent of choice, which should be purged with an inert gas. DS44960156 is soluble in organic solvents such as DMSO and methanol.

Description

DS44960156 is an inhibitor of methylenetetrahydrofolate dehydrogenase 2 (MTHFD2; IC₅₀ = 1.6 μM).¹ It is selective for MTHFD2 over MTHFD1 (IC₅₀ = >30 μM).

Reference

1. Kawai, J., Ota, M., Ohki, H., *et al.* Structure-based design and synthesis of an isozyme-selective MTHFD2 inhibitor with a tricyclic coumarin scaffold. *ACS Med. Chem. Lett.* **10(6)**, 893-898 (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, 11/09/2020

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