

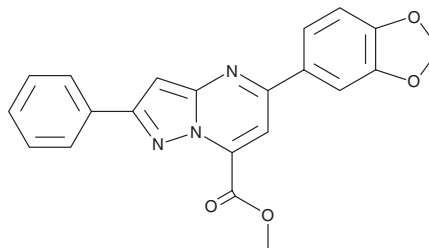
PRODUCT INFORMATION



MCI-INI-3

Item No. 42077

CAS Registry No.: 1011366-31-6
Formal Name: 5-(1,3-benzodioxol-5-yl)-2-phenyl-pyrazolo[1,5-a]pyrimidine-7-carboxylic acid, methyl ester
MF: C₂₁H₁₅N₃O₄
FW: 373.4
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

MCI-INI-3 is supplied as a solid. A stock solution may be made by dissolving the MCI-INI-3 in the solvent of choice, which should be purged with an inert gas. MCI-INI-3 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

MCI-INI-3 is an inhibitor of aldehyde dehydrogenase 1A3 (ALDH1A3; $K_i = 0.55 \mu\text{M}$).¹ It is selective for ALDH1A3 over ALDH1A1 ($K_i = 78.2 \mu\text{M}$). MCI-INI-3 (15 μM) inhibits the conversion of retinol to retinaldehyde in U87MG glioblastoma cells. It reduces the viability of GSC-83 and GSC-326 glioblastoma cells when used at a concentration of 15 μM .

Reference

1. Li, J., Garavaglia, S., Ye, Z., *et al.* A specific inhibitor of ALDH1A3 regulates retinoic acid biosynthesis in glioma stem cells. *Commun. Biol.* **4**(1), 1420 (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, 10/08/2024

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