

# PRODUCT INFORMATION

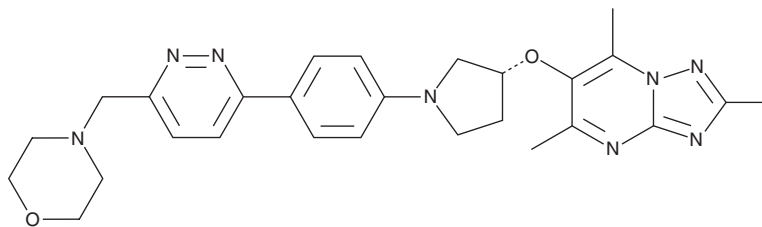


**AZD 0095**

Item No. 42134

**CAS Registry No.:** 2750001-23-9  
**Formal Name:** 2,5,7-trimethyl-6-[[[(3R)-1-[4-[6-(4-morpholinylmethyl)-3-pyridazinyl]phenyl]-3-pyrrolidinyl]oxy]-[1,2,4]triazolo[1,5-a]pyrimidine

**MF:** C<sub>27</sub>H<sub>32</sub>N<sub>8</sub>O<sub>2</sub>  
**FW:** 500.6  
**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

AZD 0095 is supplied as a solid. A stock solution may be made by dissolving the AZD 0095 in the solvent of choice, which should be purged with an inert gas. AZD 0095 is sparingly soluble (1-10 mg/ml) in DMSO.

## Description

AZD 0095 is an inhibitor of monocarboxylate transporter 4 (MCT4; IC<sub>50</sub> = 1.3 nM for lactate efflux in SK-BR-3 cells).<sup>1</sup> It selectively inhibits lactate influx in NCI H358 non-small cell lung cancer cells (IC<sub>50</sub> = 6.1 nM), which highly express MCT4 and do not express MCT1, over lactate influx in K562 chronic myeloid leukemia cells (IC<sub>50</sub> = >30 μM), which highly express MCT1 and do not express MCT4. It inhibits the proliferation of NCI H358 cells with an IC<sub>50</sub> value of 26 nM. AZD 0095 (100 mg/kg twice per day) in combination with cediranib (Item No. 11495) reduces tumor growth in an NCI H358 mouse xenograft model.

## Reference

1. Goldberg, F.W., Kettle, J.G., Lamont, G.M., *et al.* Discovery of clinical candidate AZD0095, a selective inhibitor of monocarboxylate transporter 4 (MCT4) for oncology. *J. Med. Chem.* **66**(1), 384-397 (2023).

### 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, 11/07/2024

## 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