

PRODUCT INFORMATION

Glutaminase C Inhibitor 11

Item No. 39883

CAS Registry No.: 3056437-77-2
Formal Name: 3-phenoxy-N-[5-[[[(3R)-1-(3-pyridazinyl)-3-pyrrolidinyl]amino]-1,3,4-thiadiazol-2-yl]-benzeneacetamide

MF: C₂₄H₂₃N₇O₂S

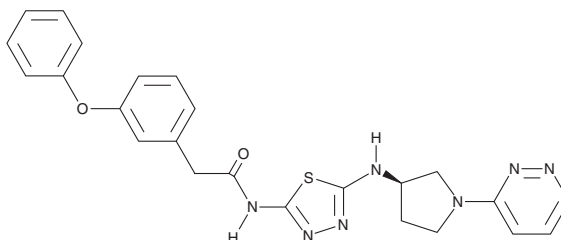
FW: 473.6

Purity: ≥95%

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

Glutaminase C inhibitor 11 is supplied as a solid. A stock solution may be made by dissolving the glutaminase C inhibitor 11 in the solvent of choice, which should be purged with an inert gas. Glutaminase C inhibitor 11 is soluble in the organic solvent DMSO.

Description

Glutaminase C inhibitor 11 is an inhibitor of glutaminase C (GAC).¹ It inhibits GAC enzyme activity (EC₅₀ = 10.64 nM) and reduces the viability of A549 non-small cell lung cancer (NSCLC) cells (IC₅₀ = 4.025 nM). Glutaminase C inhibitor 11 (0.1 and 1 μM) reduces colony formation by, and intracellular glutamate levels in, A549 cells. *In vivo*, glutaminase C inhibitor 11 (100 mg/kg) reduces tumor growth in an A549 mouse xenograft model.

Reference

1. Sun, H., Du, T., Yang, M., *et al.* Targeting the subpocket enables the discovery of thiadiazole-pyridazine derivatives as glutaminase C inhibitors. *ACS Med. Chem. Lett.* **14**(10), 1455-1466 (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, 06/19/2025

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