

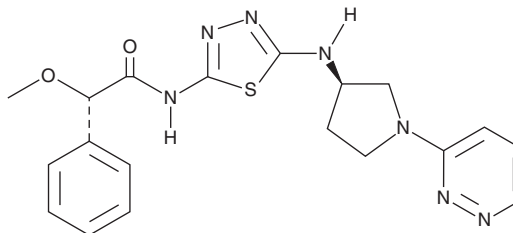
PRODUCT INFORMATION



GLS1 Inhibitor

Item No. 28581

CAS Registry No.: 1832646-52-2
Formal Name: αS-methoxy-N-[5-[[[(3R)-1-(3-pyridazinyl)-3-pyrrolidinyl]amino]-1,3,4-thiadiazol-2-yl]-benzeneacetamide
MF: C₁₉H₂₁N₇O₂S
FW: 411.5
Purity: ≥98%
UV/Vis.: λ_{max}: 252, 289 nm
Supplied as: A crystalline 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

GLS1 inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the GLS1 inhibitor in the solvent of choice. GLS1 inhibitor is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of GLS1 inhibitor in these solvents is approximately 20 mg/ml.

GLS1 inhibitor is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GLS1 inhibitor should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GLS1 inhibitor has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GLS1 inhibitor is an inhibitor of glutaminase 1 (GLS1; IC₅₀ = 0.021 μM).¹ It inhibits the growth of NCI H1703 non-small cell lung cancer (NSCLC) cells *in vitro* (GI₅₀ = 0.011 μM). GLS1 inhibitor (100 mg/kg) reduces tumor growth, increases tumor levels of glutamine, and decreases tumor levels of glutamate and aspartate in an NCI H1703 mouse xenograft model.

Reference

1. Finlay, M.R.V., Anderton, M., Bailey, A., *et al.* Discovery of a thiadiazole-pyridazine-based allosteric glutaminase 1 inhibitor series that demonstrates oral bioavailability and activity in tumor xenograft models. *J. Med. Chem.* **62**(14), 6540-6560 (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, 03/19/2024

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