

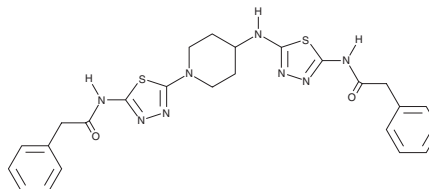
# PRODUCT INFORMATION



**UPGL00004**

Item No. 42323

**CAS Registry No.:** 1890169-95-5  
**Formal Name:** N-[5-[4-[[5-[(2-phenylacetyl)amino]-1,3,4-thiadiazol-2-yl]amino]-1-piperidiny]-1,3,4-thiadiazol-2-yl]-benzeneacetamide  
**Synonym:** UP4  
**MF:** C<sub>25</sub>H<sub>26</sub>N<sub>8</sub>O<sub>2</sub>S<sub>2</sub>  
**FW:** 534.7  
**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

UPGL00004 is supplied as a solid. A stock solution may be made by dissolving the UPGL00004 in the solvent of choice, which should be purged with an inert gas. UPGL00004 is soluble (≥10 mg/ml) in DMSO.

## Description

UPGL00004 is an inhibitor of glutaminase C (GAC; IC<sub>50</sub> = 29 nM).<sup>1</sup> It reduces the proliferation of MDA-MB-231, Hs 578T, and TSE breast cancer cells (IC<sub>50</sub>s = 70, 129, and 262 nM, respectively). UPGL00004 (0.1 μM) also inhibits replication of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) and human coronavirus OC43 (HCoV-OC43) in Vero and primary human bronchial epithelial cells (HBECs), respectively.<sup>2</sup>

## References

1. Huang, Q., Stalnecker, C., Zhang, C., *et al.* Characterization of the interactions of potent allosteric inhibitors with glutaminase C, a key enzyme in cancer cell glutamine metabolism. *J. Biol. Chem.* **293**(10), 3535-3545 (2018).
2. Greene, K.S., Choi, A., Yang, N., *et al.* Glutamine metabolism is essential for coronavirus replication in host cells and in mice. *J. Biol. Chem.* **301**(1), 108063 (2025).

### 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, 05/12/2025

## 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](http://WWW.CAYMANCHEM.COM)