

PRODUCT INFORMATION



Glutaminase-IN-3

Item No. 42325

CAS Registry No.: 1439399-45-7
Formal Name: N-[6-[4-(5-amino-1,3,4-thiadiazol-2-yl)butyl]-3-pyridazinyl]-3-(trifluoromethoxy)benzeneacetamide

Synonym: Glutaminase Inhibitor 3

MF: C₁₉H₁₉F₃N₆O₂S

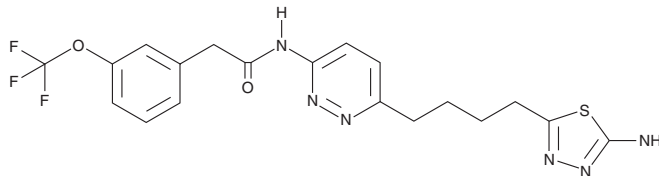
FW: 452.5

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

Glutaminase-IN-3 is supplied as a solid. A stock solution may be made by dissolving the glutaminase-IN-3 in the solvent of choice, which should be purged with an inert gas. Glutaminase-IN-3 is soluble (≥10 mg/ml) in DMSO.

Description

Glutaminase-IN-3 is an inhibitor of the glutaminase 1 (GLS1) isozyme glutaminase C (GAC; IC₅₀ = 0.24 μM).¹ It inhibits the proliferation of PC3 and LNCaP prostate cancer cells (IC₅₀s = 6.14 and 2.13 μM, respectively) and non-cancerous CCD-1072Sk fibroblasts (IC₅₀ = 15.39 μM).²

References

1. Parlati, F., Rodriguez, M.L., Gross, M.I., *et al.* Combination therapy with glutaminase inhibitors. (2015).
2. Demir, U. and Celik, A.B. The anti-proliferative effect of a novel glutaminase inhibitor IN-3 on prostate cancer cells. *Medeni Med. J.* **39(3)**, 169-174 (2024).

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, 11/06/2024

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