

PRODUCT INFORMATION



GLUT Inhibitor-1 (formate)

Item No. 42732

Formal Name: 2-[3-[6-cyclobutyl-5,6,7,8-tetrahydro-4-[[4-(1H-pyrazol-4-yl)phenyl]amino]pyrido[4,3-d]pyrimidin-2-yl]phenoxy]-N-(1-methylcyclopropyl)-acetamide, monoformic acid

Synonym: Glucose Transporter Inhibitor-1

MF: C₃₂H₃₅N₇O₂ • CH₂O₂

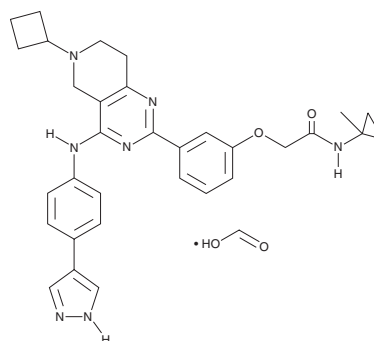
FW: 595.7

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

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

Description

GLUT inhibitor-1 is an inhibitor of glucose transporter 1 (Glut1) and Glut3 (IC₅₀s = 0.242 and 0.179 μM, respectively).¹ It is selective for Glut1 and Glut3 over a panel of 31 receptors, ion channels, and transporters at 10 μM.

Reference

1. Liu, K.G., Kim, J.-I., Olszewski, K., *et al.* Discovery and optimization of glucose uptake inhibitors. *J. Med. Chem.* **63**(10), 5201-5211 (2020).

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, 06/20/2025

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