

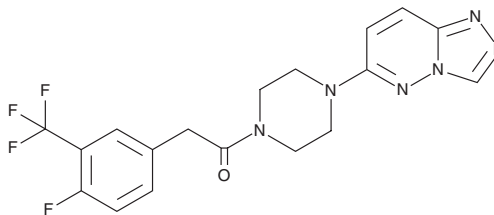
PRODUCT INFORMATION



SLU-10482

Item No. 39375

CAS Registry No.: 2755451-45-5
Formal Name: 2-[4-fluoro-3-(trifluoromethyl)phenyl]-1-[4-(1,2,4-triazolo[4,3-b]pyridazin-6-yl)-1-piperazinyl]-ethanone
MF: C₁₈H₁₆F₄N₆O
FW: 408.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SLU-10482 is supplied as a solid. A stock solution may be made by dissolving the SLU-10482 in the solvent of choice, which should be purged with an inert gas. SLU-10482 is slightly soluble in ethanol, DMSO, and methanol.

Description

SLU-10482 is an antiparasitic agent.¹ It reduces the number of *C. parvum* parasites in HCT-8 cells (EC₅₀ = 0.07 μM). SLU-10482 binds to human ether-a-go-go (hERG; K_d = 43 μM) to a lesser extent than the antiparasitic SLU-2633. *In vivo*, SLU-10482 (5 and 15 mg/kg twice per day) decreases the number of fecal oocysts in a mouse model of *C. parvum* infection.

Reference

- Schubert, T.J., Oboh, E., Peek, H., *et al.* Structure-activity relationship studies of the aryl acetamide triazolopyridazines against *Cryptosporidium* reveals remarkable role of fluorine. *J. Med. Chem.* **66**(12), 7834-7848 (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, 10/31/2023

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