

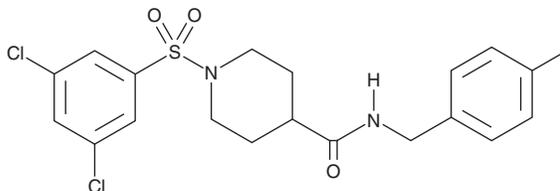
PRODUCT INFORMATION



BI-01383298

Item No. 24482

CAS Registry No.: 2227549-00-8
Formal Name: 1-[(3,5-dichlorophenyl)sulfonyl]-
N-[(4-fluorophenyl)methyl]-4-
piperidinecarboxamide
MF: C₁₉H₁₉Cl₂FN₂O₃S
FW: 445.3
Purity: ≥95%
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

BI-01383298 is supplied as a crystalline solid. A stock solution may be made by dissolving the BI-01383298 in the solvent of choice, which should be purged with an inert gas. BI-01383298 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BI-01383298 in these solvents is approximately 5 and 30 mg/ml, respectively.

Description

BI-01383298 is an inhibitor of the sodium-coupled citrate transporter (NaCT; IC₅₀ = ~100 nM).¹ It irreversibly inhibits sodium-dependent citrate uptake, but not nicotinate, arginine, glutamine, tryptophan, or lactate uptake, in HepG2 cells when used at concentrations ranging from 0.01 to 1 μM.

Reference

1. Higuchi, K., Kopel, J.J., Sivaprakasam, S., *et al.* Functional analysis of a species-specific inhibitor selective for human Na⁺-coupled citrate transporter (NaCT/SLC13A5/mINDY). *Biochem. J.* **477(21)**, 4149-4165 (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, 11/22/2022

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