

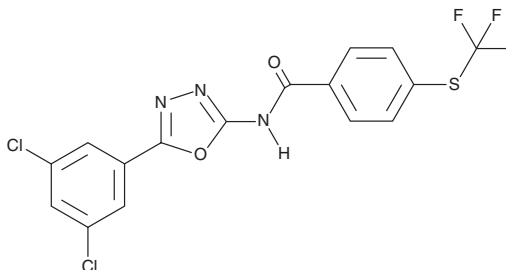
PRODUCT INFORMATION



HSGN-218

Item No. 32711

CAS Registry No.: 2519410-44-5
Formal Name: N-[5-(3,5-dichlorophenyl)-1,3,4-oxadiazol-2-yl]-4-[(trifluoromethyl)thio]-benzamide
MF: C₁₆H₈Cl₂F₃N₃O₂S
FW: 434.2
Purity: ≥98%
UV/Vis.: λ_{max}: 281 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

HSGN-218 is supplied as a crystalline solid. A stock solution may be made by dissolving the HSGN-218 in the solvent of choice, which should be purged with an inert gas. HSGN-218 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HSGN-218 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

HSGN-218 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HSGN-218 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. HSGN-218 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HSGN-218 is a gut-restrictive antibiotic.¹ It is active against *C. difficile* (MIC = 0.007 µg/ml) and clinical isolates of *C. difficile* (MICs = 0.003-0.03 µg/ml) and is not cytotoxic to Caco-2 cells when used at concentrations ranging from 4 to 64 µg/ml. HSGN-218 (50 mg/kg) increases survival and protects mice from *C. difficile* recurrence in mouse models of *C. difficile* infection.

Reference

1. Naclerio, G.A., Abutaleb, N.S., Li, D., *et al.* Ultrapotent inhibitor of *Clostridioides difficile* growth, which suppresses recurrence *in vivo*. *J. Med. Chem.* **63**(20), 11934-11944 (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, 07/27/2023

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