

PRODUCT INFORMATION



Voxilaprevir

Item No. 33974

CAS Registry No.: 1535212-07-7
Formal Name: N-[[[(1R,2R)-2-[5,5-difluoro-5-(3-hydroxy-6-methoxy-2-quinoxaliny)pentyl]cyclopropyl]oxy]carbonyl]-3-methyl-L-valyl-(3S,4R)-3-ethyl-4-hydroxy-L-prolyl-1-amino-2-(difluoromethyl)-N-[(1-methylcyclopropyl)sulfonyl]-cyclopropanecarboxamide, cyclic (1R→2R)-ether GS-9857

Synonym:

MF: C₄₀H₅₂F₄N₆O₉S

FW: 868.9

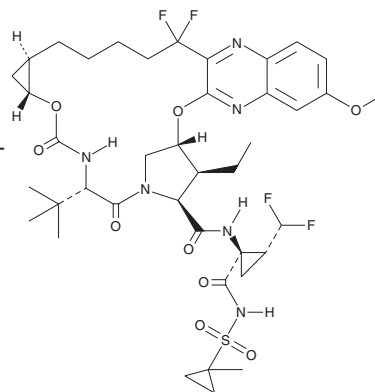
Purity: ≥98%

UV/Vis.: λ_{max}: 217 nm

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

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

Description

Voxilaprevir is a pan-genotypic hepatitis C virus (HCV) non-structural 3/4A (NS3/4A) serine protease inhibitor.¹ It inhibits replication of wild-type HCV genotypes 1a, 1b, 2a, 2b, 3a, 4a, 5a, and 6a (EC₅₀s = 1.9-6.6 nM in replicon assays). Voxilaprevir also inhibits replication of mutant HCV genotypes containing NS3 resistance-associated substitutions (EC₅₀s = 3.1, 2.7, and 4 nM for HCV genotypes 1a Q80K, 1a R155K, and 1b D168E, respectively). Formulations containing voxilaprevir have been used in the treatment of chronic HCV infection.

Reference

1. Taylor, J.G., Zipfel, S., Ramey, K., *et al.* Discovery of the pan-genotypic hepatitis C virus NS3/4A protease inhibitor voxilaprevir (GS-9857): A component of Vosevi®. *Bioorg. Med. Chem. Lett.* **29(16)**, 2428-2436 (2019).

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, 10/10/2022

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