

PRODUCT INFORMATION

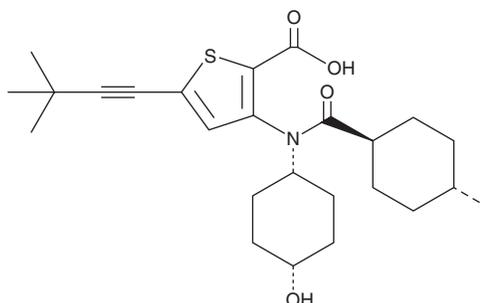


Lomibuvir

Item No. 26630

CAS Registry No.: 1026785-59-0
Formal Name: 5-(3,3-dimethyl-1-butyne-1-yl)-3-[[*cis*-4-hydroxycyclohexyl][[*trans*-4-methylcyclohexyl]carbonyl]amino]-2-thiophenecarboxylic acid

Synonym: VX-222
MF: C₂₅H₃₅NO₄S
FW: 445.6
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 299 nm
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

Lomibuvir is supplied as a crystalline solid. A stock solution may be made by dissolving the lomibuvir in the solvent of choice, which should be purged with an inert gas. Lomibuvir is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

Lomibuvir is a non-nucleoside inhibitor of the RNA-directed RNA polymerase of hepatitis C virus (HCV; K_i = 17 nM) that inhibits primer-dependent RNA synthesis selectively over *de novo* RNA synthesis.¹ It inhibits HCV replication in Huh7.5 cells expressing the 1b/Con1 HCV subgenomic replicon (EC₅₀ = 5 nM).

Reference

1. Yi, G., Deval, J., Fan, B., *et al.* Biochemical study of the comparative inhibition of hepatitis C virus RNA polymerase by VX-222 and filibuvir. *Antimicrob. Agents Chemother.* **56(2)**, 830-837 (2012).

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, 11/02/2022

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