

PRODUCT INFORMATION



Coblopasvir (hydrochloride)

Item No. 38793

CAS Registry No.: 1966138-53-3
Formal Name: N-[(1S)-1-[[[(2S)-2-[5-[4-[7-[2-[(2S)-1-[(2S)-2-[(methoxycarbonyl)amino]-3-methyl-1-oxobutyl]-2-pyrrolidinyl]-1H-imidazol-5-yl]-1,3-benzodioxol-4-yl]phenyl]-1H-imidazol-2-yl]-1-pyrrolidinyl]carbonyl]-2-methylpropyl]-carbamic acid, methyl ester, dihydrochloride

MF: C₄₁H₅₀N₈O₈ • 2HCl

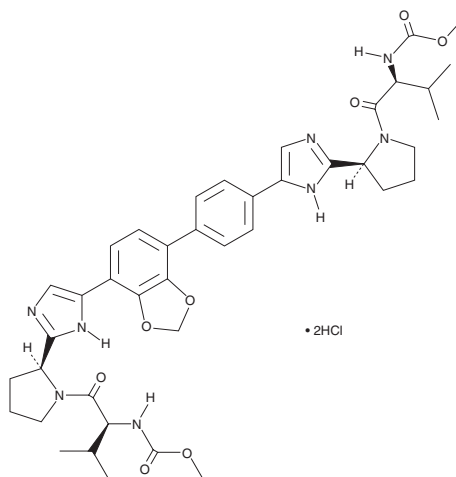
FW: 855.8

Purity: ≥98%

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

Coblopasvir (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the coblopasvir (hydrochloride) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Coblopasvir is a pan-genotypic inhibitor of hepatitis C virus (HCV) non-structural protein 5A (NS5A).¹ It inhibits viral replication of the HCV genotypes 1a, -1b, -3a, -4a, -5a, and -6a in cellular assays and genotype 2a in a cell-free assay.

Reference

1. Rao, H., Song, G., Li, G., *et al.* Safety and efficacy of coblopasvir and sofosbuvir in patients with genotypes 1, 2, 3 and 6 HCV infections without or with compensated cirrhosis. *J. Viral. Hepat.* **27**(1), 45-51 (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, 07/26/2023

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