

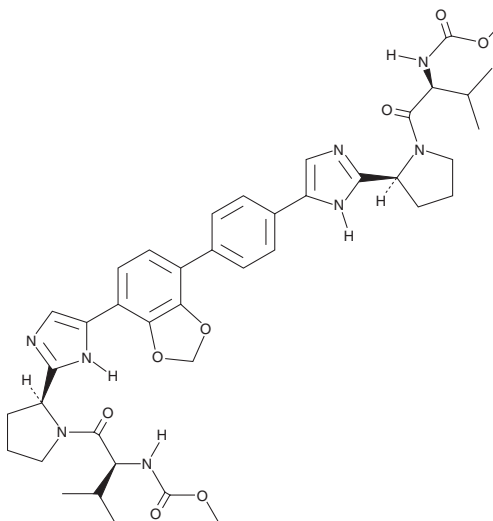
# PRODUCT INFORMATION



## Coblopassvir

Item No. 36779

**CAS Registry No.:** 1312608-46-0  
**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  
**Synonym:** KW-136  
**MF:** C<sub>41</sub>H<sub>50</sub>N<sub>8</sub>O<sub>8</sub>  
**FW:** 782.9  
**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

Coblopassvir is supplied as a solid. A stock solution may be made by dissolving the coblopassvir in the solvent of choice, which should be purged with an inert gas. Coblopassvir is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of coblopassvir in these solvents is approximately 10 and 3 mg/ml, respectively. Coblopassvir is also slightly soluble in ethanol.

### Description

Coblopassvir is a pan-genotypic inhibitor of hepatitis C virus (HCV) non-structural protein 5A (NS5A).<sup>1</sup> 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 coblopassvir 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, 12/21/2022

#### 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](http://WWW.CAYMANCHEM.COM)