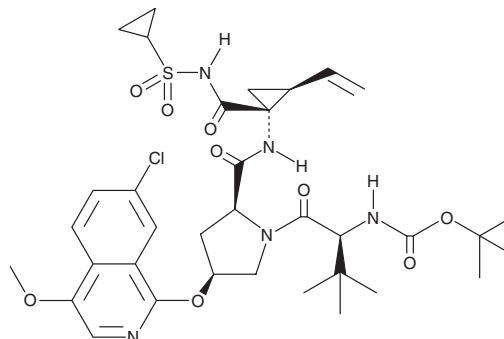


# PRODUCT INFORMATION



## Asunaprevir Item No. 20835

**CAS Registry No.:** 630420-16-5  
**Formal Name:** (1R,2S)-N-[(1,1-dimethylethoxy)carbonyl]-3-methyl-L-valyl-(4R)-4-[(7-chloro-4-methoxy-1-isoquinolinyl)oxy]-L-prolyl-1-amino-N-(cyclopropylsulfonyl)-2-ethenyl-cyclopropanecarboxamide  
**Synonym:** BMS 650032  
**MF:** C<sub>35</sub>H<sub>46</sub>ClN<sub>5</sub>O<sub>9</sub>S  
**FW:** 748.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 254, 305, 333 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

Asunaprevir is supplied as a crystalline solid. A stock solution may be made by dissolving the asunaprevir in the solvent of choice, which should be purged with an inert gas. Asunaprevir is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of asunaprevir in these solvents is approximately 15, 25, and 30 mg/ml, respectively.

Asunaprevir is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, asunaprevir should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Asunaprevir has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Asunaprevir is an inhibitor of the hepatitis C non-structural 3/4A serine protease, which is required for viral replication.<sup>1</sup> It is a direct-acting inhibitor of hepatitis C virus (HCV).<sup>1,2</sup> It is commonly used in combination with daclatasvir and beclabuvir, which are also direct-acting HCV inhibitors.<sup>1</sup>

### References

1. Zhang, X. Direct anti-HCV agents *Acta Pharmaceutica Sinica B* **6(1)**, 26-31 (2015).
2. Alavian, S.M. and Rezaee-Zavareh, M.S. Daclatasvir-based treatment regimens for hepatitis C virus infection: A systematic review and meta-analysis. *Hepat. Mon.* **16(9)**, e41077 (2016).

#### 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, 11/09/2022

#### 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