

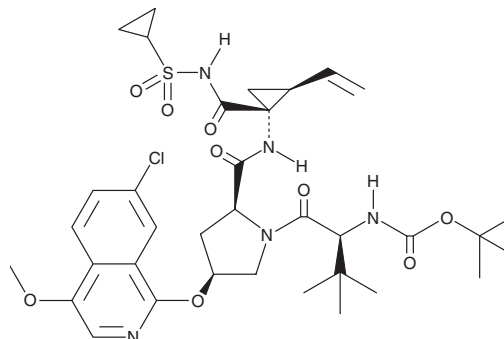
PRODUCT INFORMATION



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-ethenylcyclopropanecarboxamide

Synonym: BMS 650032
MF: C₃₅H₄₆ClN₅O₉S
FW: 748.3
Purity: ≥98%
UV/Vis.: λ_{max}: 254, 305, 333 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

Asunaprevir is supplied as a crystalline solid. A stock solution may be made by dissolving the asunaprevir in the solvent of choice. Asunaprevir is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. 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.¹ It is a direct-acting inhibitor of hepatitis C virus (HCV).^{1,2} It is commonly used in combination with daclatasvir and beclabuvir, which are also direct-acting HCV inhibitors.¹

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, 01/03/2017

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