

PRODUCT INFORMATION

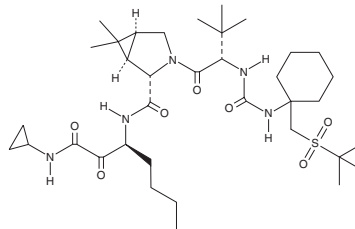


Narlaprevir

Item No. 42487

CAS Registry No.: 865466-24-6
Formal Name: (1R,2S,5S)-N-[(1S)-1-[2-(cyclopropylamino)-2-oxoacetyl]pentyl]-3-[(2S)-2-[[[1-[[[(1,1-dimethylethyl)sulfonyl]methyl]cyclohexyl]amino]carbonyl]amino]-3,3-dimethyl-1-oxobutyl]-6,6-dimethyl-3-azabicyclo[3.1.0]hexane-2-carboxamide

Synonym: SCH 900518
MF: C₃₆H₆₁N₅O₇S
FW: 708.0
Purity: ≥95%
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

Narlaprevir is supplied as a solid. A stock solution may be made by dissolving the narlaprevir in the solvent of choice, which should be purged with an inert gas. Narlaprevir is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Narlaprevir is an inhibitor of hepatitis C virus (HCV) non-structural protein 3 (NS3; $K_i = 6$ nM).¹ It also inhibits severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) main protease (M^{pro}), also known as 3C-like protease (3CL^{pro}; $IC_{50} = 16.11$ μM).² Narlaprevir inhibits HCV replication in a cell-based replicon assay ($EC_{90} = 40$ nM) and SARS-CoV-2 replication in Vero cells ($EC_{50} = 7.43$ μM).^{1,2}

References

1. Arasappan, A., Bennett, F., Bogen, S.L., *et al.* Discovery of narlaprevir (SCH 900518): A potent, second generation HCV NS3 serine protease inhibitor. *ACS Med. Chem. Lett.* **1(2)**, 64-69 (2010).
2. Bai, Y., Ye, F., Feng, Y., *et al.* Structural basis for the inhibition of the SARS-CoV-2 main protease by the anti-HCV drug narlaprevir. *Signal Transduct. Target. Ther.* **6(1)**, 51 (2024).

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, 12/18/2024

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