

PRODUCT INFORMATION



Samatasvir

Item No. 36776

CAS Registry No.: 1312547-19-5
Formal Name: N-[(1R)-2-[(2S)-2-[5-[4-[6-[2-[(2S)-1-[(2S)-2-[(methoxycarbonyl)amino]-3-methyl-1-oxobutyl]-2-pyrrolidinyl]-1H-benzimidazol-6-yl]thieno[3,2-b]thien-3-yl]phenyl]-1H-imidazol-2-yl]-1-pyrrolidinyl]-2-oxo-1-phenylethyl]-carbamic acid, methyl ester

Synonyms: IDX719, IDX18719

MF: C₄₇H₄₈N₈O₆S₂

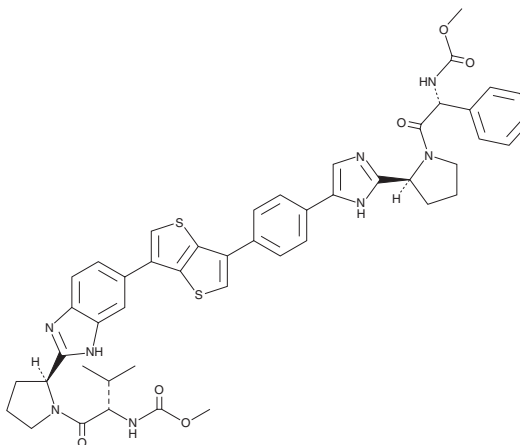
FW: 885.1

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

Samatasvir is supplied as a solid. A stock solution may be made by dissolving the samatasvir in the solvent of choice, which should be purged with an inert gas. Samatasvir is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of samatasvir in these solvents is approximately 1 and 3 mg/ml, respectively.

Description

Samatasvir is an inhibitor of hepatitis C virus (HCV) non-structural protein 5A (NS5A).¹ It inhibits HCV replication in stable replicon cell lines containing NS5A from genotypes 1a, -1b, -2a, -3a, -4a, and -5a (EC₅₀s = 2-24 pM). Samatasvir acts additively with the nucleotide prodrug IDX184, NS5B inhibitor TMC647055, or NS3/4A protease inhibitor simeprevir (Item No. 22144) to inhibit viral replication in genotype 1b HCV replicon cells.

Reference

1. Bilello, J.P., Lалlos, L.B., McCarville, J.F., *et al.* *In vitro* activity and resistance profile of samatasvir, a novel NS5A replication inhibitor of hepatitis C virus. *Antimicrob. Agents Chemother.* **58(8)**, 4431-4442 (2014).

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