

PRODUCT INFORMATION



Ravidasvir (hydrochloride)

Item No. 37794

CAS Registry No.: 1303533-81-4

Formal Name: N-[(1S)-1-[[[(2S)-2-[5-[6-[2-[(2S)-1-[(2S)-2-[(methoxycarbonyl)amino]-3-methyl-1-oxobutyl]-2-pyrrolidinyl]-1H-benzimidazol-6-yl]-2-naphthalenyl]-1H-imidazol-2-yl]-1-pyrrolidinyl]carbonyl]-2-methylpropyl]-carbamic acid, methyl ester, dihydrochloride

Synonym: PPI-668

MF: $C_{42}H_{50}N_8O_6 \cdot 2HCl$

FW: 835.8

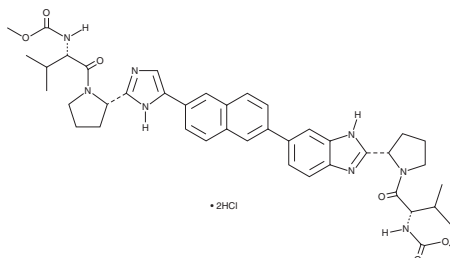
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 267, 321 nm

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ravidasvir (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the ravidasvir (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Ravidasvir (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ravidasvir (hydrochloride) in ethanol is approximately 1 mg/ml and approximately 5 mg/ml in DMSO and DMF.

Description

Ravidasvir is an inhibitor of hepatitis C virus (HCV) non-structural protein 5A (NS5A).¹ It inhibits HCV replication in Huh7 replicon cells containing NS5A from the HCV genotypes 1a, -1b, and 3a with EC_{50} values of 0.12, 0.01, and 1.14 nM, respectively.

Reference

1. Zhong, M., Peng, E., Huang, N., *et al.* Discovery of ravidasvir (PPI-668) as a potent pan-genotypic HCV NS5A inhibitor. *Bioorg. Med. Chem. Lett.* **26(18)**, 4508-4512 (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, 06/05/2023

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