

PRODUCT INFORMATION

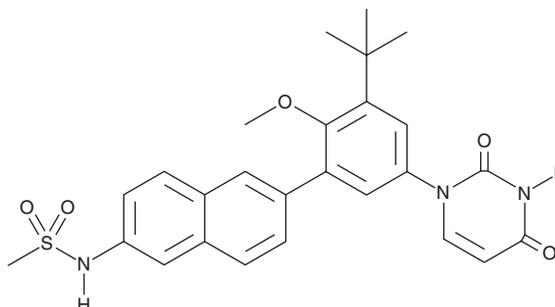


ABT-333

Item No. 18482

CAS Registry No.: 1132935-63-7
Formal Name: N-[6-[5-(3,4-dihydro-2,4-dioxo-1(2H)-pyrimidinyl)-3-(1,1-dimethylethyl)-2-methoxyphenyl]-2-naphthalenyl]-methanesulfonamide

Synonym: Dasabuvir
MF: C₂₆H₂₇N₃O₅S
FW: 493.5
Purity: ≥98%
UV/Vis.: λ_{max}: 244 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

ABT-333 is supplied as a crystalline solid. A stock solution may be made by dissolving the ABT-333 in the solvent of choice, which should be purged with an inert gas. ABT-333 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of ABT-333 in these solvents is approximately 20 and 30 mg/ml, respectively.

ABT-333 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ABT-333 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ABT-333 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ABT-333 is an inhibitor of hepatitis C virus (HCV) RNA-dependent RNA polymerase (RdRp; IC₅₀s = 0.0022-0.0107 μM for genotypes 1a and 1b).¹ It is selective for HCV genotype 1 polymerases over genotypes 2, 3, and 4 (IC₅₀s = 0.9->20 μM), as well as human polymerases (IC₅₀s = >100 μM). Formulations containing ABT-333, ombitasvir, paritaprevir, and ritonavir have been used in the treatment of HCV genotype 1 infections.

Reference

1. Kati, W., Koev, G., Irvin, M., *et al.* *In vitro* activity and resistance profile of dasabuvir, a nonnucleoside hepatitis C virus polymerase inhibitor. *Antimicrob. Agents Chemother.* **59**(3), 1505-1511 (2015).

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