

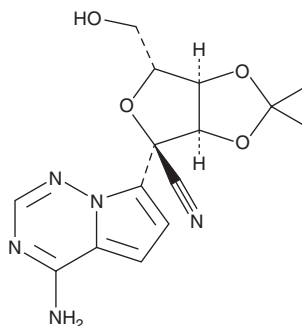
PRODUCT INFORMATION



Remdesivir O-desphosphate acetonide

Item No. 36673

CAS Registry No.:	1191237-80-5
Formal Name:	2-C-(4-aminopyrrolo[2,1-f][1,2,4]triazin-7-yl)-2,5-anhydro-3,4-O-(1-methylethylidene)-D-altronitrile
MF:	C ₁₅ H ₁₇ N ₅ O ₄
FW:	331.3
Purity:	≥98%
UV/Vis.:	λ _{max} : 245 nm
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

Remdesivir O-desphosphate acetonide is supplied as a solid. A stock solution may be made by dissolving the remdesivir O-desphosphate acetonide in the solvent of choice, which should be purged with an inert gas. Remdesivir O-desphosphate acetonide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of remdesivir O-desphosphate acetonide in these solvents is approximately 5, 11, and 12 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of remdesivir O-desphosphate acetonide can be prepared by directly dissolving the solid in aqueous buffers. The solubility of remdesivir O-desphosphate acetonide in PBS (pH 7.2) is approximately 0.30 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Remdesivir O-desphosphate acetonide is a synthetic intermediate in the synthesis of the antiviral prodrug remdesivir (Item No. 30354).¹

Reference

1. Gannedi, V., Villuri, B.K., Reddy, S.N., *et al.* Practical remdesivir synthesis through one-pot organocatalyzed asymmetric (S)-P-phosphoramidation. *J. Org. Chem.* **86(7)**, 4977-4985 (2021).

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, 10/05/2022

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