

# PRODUCT INFORMATION



## Remdesivir isopropyl ester analog

Item No. 9003635

**CAS Registry No.:** 1911578-83-0  
**Formal Name:** N-[(S)-hydroxyphenoxyphosphinyl]-L-alanine, 1-methylethyl ester, 6-ester with 2-C-(4-aminopyrrolo[2,1-f][1,2,4]triazin-7-yl)-2,5-anhydro-D-altrnonitrile

**Synonyms:** Anti-virus agent 1,  
GS-5734 isopropyl ester analog

**MF:** C<sub>24</sub>H<sub>29</sub>N<sub>6</sub>O<sub>8</sub>P

**FW:** 560.5

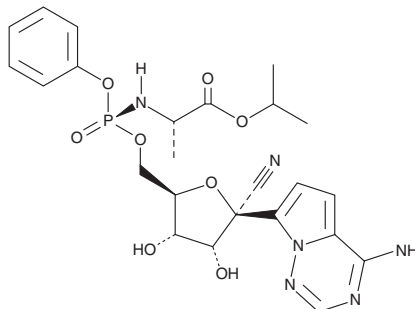
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 246 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

Remdesivir isopropyl ester analog is supplied as a crystalline solid. A stock solution may be made by dissolving the remdesivir isopropyl ester analog in the solvent of choice, which should be purged with an inert gas. Remdesivir isopropyl ester analog is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of remdesivir isopropyl ester analog in these solvents is approximately 10 mg/ml.

Remdesivir isopropyl ester analog is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, remdesivir isopropyl ester analog should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Remdesivir isopropyl ester analog has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Remdesivir isopropyl ester analog is a derivative of the antiviral prodrug remdesivir (Item No. 30354).

#### 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, 05/02/2021

#### 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](http://WWW.CAYMANCHEM.COM)