

# PRODUCT INFORMATION



## Remdesivir methylpropyl ester analog

Item No. 9003571

**Formal Name:** N-[(S)-hydroxyphenoxyphosphinyl]-L-alanine, 2-methylpropyl ester, 6-ester with 2-C-(4-aminopyrrolo[2,1-f][1,2,4]triazin-7-yl)-2,5-anhydro-D-altronitrile

**Synonym:** GS-5734 methylpropyl ester analog

**MF:** C<sub>25</sub>H<sub>31</sub>N<sub>6</sub>O<sub>8</sub>P

**FW:** 574.5

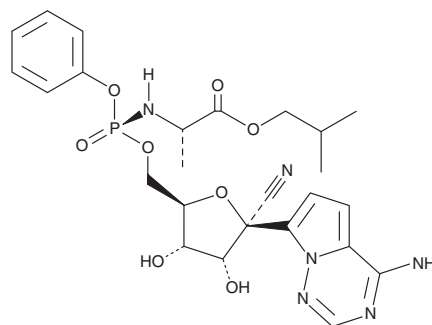
**Purity:** ≥98%

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

Remdesivir methylpropyl ester analog is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, remdesivir methylpropyl ester analog should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Remdesivir methylpropyl 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 methylpropyl 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, 05/02/2024

#### 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)