

# PRODUCT INFORMATION



**GS-704277**

Item No. 9003569

**CAS Registry No.:** 1911579-04-8

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

**MF:** C<sub>15</sub>H<sub>19</sub>N<sub>6</sub>O<sub>8</sub>P

**FW:** 442.3

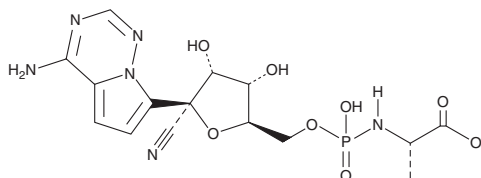
**Purity:** ≥95%

**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

GS-704277 is supplied as a crystalline solid. A stock solution may be made by dissolving the GS-704277 in the solvent of choice, which should be purged with an inert gas. GS-704277 is soluble in the organic solvent DMSO.

## Description

GS-704277 is an intermediate alanine metabolite of the antiviral prodrug remdesivir (Item No. 30354).<sup>1,2</sup> It is formed from remdesivir *via* intracellular hydrolysis. GS-704277 is further metabolized to the nucleoside monophosphate form and then to the nucleoside analog GS-441524 (Item No. 30469) or active nucleoside triphosphate GS-443902.

## References

1. Humeniuk, R., Mathias, A., Cao, H., *et al.* Safety, tolerability, and pharmacokinetics of remdesivir, an antiviral for treatment of COVID-19, in healthy subjects. *Clin. Transl. Sci.* **13**(5), 896-906 (2020).
2. Eastman, R.T., Roth, J.S., Brimacombe, K.R., *et al.* Remdesivir: A review of its discovery and development leading to emergency use authorization for treatment of COVID-19. *ACS Cent. Sci.* **6**(5), 672-683 (2020).

### 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, 05/04/2024

## 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