

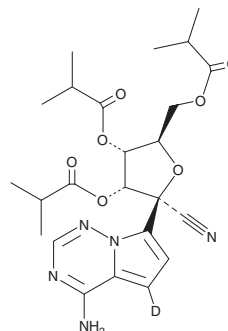
PRODUCT INFORMATION



VV116

Item No. 37745

CAS Registry No.: 2647442-33-7
Formal Name: (2R,3R,4R,5R)-2-(4-aminopyrrolo[2,1-f][1,2,4]triazin-7-yl-5-d)-2-cyano-5-((isobutyryloxy)methyl)tetrahydrofuran-3,4-diyl, bis(2-methylpropanoate)
Synonym: JT001
MF: C₂₄H₃₀DN₅O₇
FW: 502.5
Purity: ≥95%
UV/Vis.: λ_{max}: 244 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

VV116 is supplied as a solid. A stock solution may be made by dissolving the VV116 in the solvent of choice, which should be purged with an inert gas. VV116 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of VV116 in these solvents is approximately 12 and 20 mg/ml, respectively. VV116 is slightly soluble in ethanol.

Description

VV116 is an antiviral agent and a derivative of the antiviral prodrug remdesivir (Item Nos. 30354 | 32511).¹ It inhibits viral replication of respiratory syncytial virus (RSV) in A549 cells (EC₅₀ = 1.01 μM). VV116 (25 mg/kg) decreases viral replication and lung viral titer counts in a mouse model of RSV infection. It also decreases alveolar inflammation in RSV-infected mice.

Reference

1. Zhang, R., Zhang, Y., Zheng, W., *et al.* Oral remdesivir derivative VV116 is a potent inhibitor of respiratory syncytial virus with efficacy in mouse model. *Signal Transduct. Target. Ther.* **7**(1), 123 (2022).

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