

PRODUCT INFORMATION

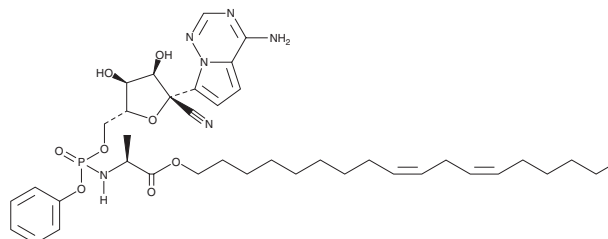


(±)-MMT5-14

Item No. 37829

CAS Registry No.: 2719679-31-7
Formal Name: (9Z,12Z)-octadeca-9,12-dien-1-yl (((2R,3S,4R,5R)-5-(4-aminopyrrolo[2,1-f][1,2,4]triazin-7-yl)-5-cyano-3,4-dihydroxytetrahydrofuran-2-yl)methoxy)(phenoxy)phosphoryl)-L-alaninate

MF: C₃₉H₅₅N₆O₈P
FW: 766.9
Purity: ≥98%
UV/Vis.: λ_{max}: 242, 246 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

(±)-MMT5-14 is supplied as a solid. A stock solution may be made by dissolving the (±)-MMT5-14 in the solvent of choice, which should be purged with an inert gas. (±)-MMT5-14 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (±)-MMT5-14 in DMF is approximately 20 mg/ml and approximately 16 mg/ml in ethanol and DMSO.

Description

(±)-MMT5-14 is a derivative of remdesivir (Item Nos. 32511 | 30354) with antiviral activity.¹ It increases the viability of Vero E6 cells infected with severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2; EC₅₀ = 0.4 μM), as well as Vero E6 cells infected with the Alpha, Beta, Gamma, or Delta SARS-CoV-2 variants (EC₅₀s = 2.5, 15.9, 1.7, and 5.6 μM, respectively). (±)-MMT5-14 exhibits increased plasma and lung concentrations compared with remdesivir in hamsters when administered at a dose of 10 mg/kg.

Reference

1. Hu, H., Traore, M.D.M., Li, R., *et al.* Optimization of the prodrug moiety of remdesivir to improve lung exposure/selectivity and enhance anti-SARS-CoV-2 activity. *J. Med. Chem.* **65**(18), 12044-12054 (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, 05/08/2024

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