

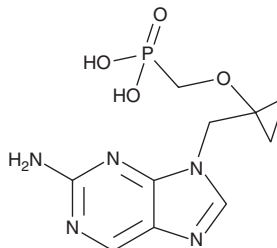
PRODUCT INFORMATION



Besifovir

Item No. 37570

CAS Registry No.: 441785-25-7
Formal Name: P-[[[1-[(2-amino-9H-purin-9-yl)methyl]cyclopropyl]oxy]methyl]-phosphonic acid
Synonym: BFV
MF: C₁₀H₁₄N₅O₄P
FW: 299.2
Purity: ≥95%
UV/Vis.: λ_{max}: 226 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

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

Description

Besifovir is an antiviral nucleoside derivative of 2'-deoxyguanosine 5'-monophosphate (dGMP; Item No. 16408).¹ It reduces viral replication in Huh7 cells infected with wild-type hepatitis B virus (HBV) or a lamivudine-resistant HBV clinical isolate (IC₅₀s = 3.5 and 7.47 μM, respectively). Besifovir also reduces viral replication in Huh7 cells infected with an entecavir-resistant HBV clinical isolate (IC₅₀ = 26 μM).

Reference

1. Won, J., Lee, A.R., Dezhbord, M., *et al.* Susceptibility of drug resistant hepatitis B virus mutants to besifovir. *Biomedicines* **10**(7), 1637 (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, 05/15/2023

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