

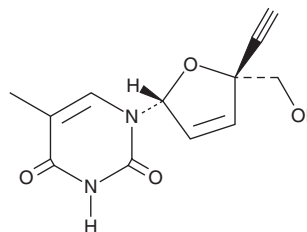
PRODUCT INFORMATION



Festinavir

Item No. 37668

CAS Registry No.: 634907-30-5
Formal Name: 2',3'-didehydro-3'-deoxy-4'-ethynyl-thymidine
Synonyms: BMS 986001, 4'-Ed4T
MF: C₁₂H₁₂N₂O₄
FW: 248.2
Purity: ≥95%
UV/Vis.: λ_{max}: 210, 265 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

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

Description

Festinavir is a nucleoside reverse transcriptase inhibitor (NRTI).¹ It inhibits viral replication in HEK293 cells infected with HIV-1 expressing wild-type or M184V mutant reverse transcriptase (EC₅₀s = 0.323 and 2.3 μM, respectively).

Reference

1. Li, Z., Terry, B., Olds, W., *et al.* *In vitro* cross-resistance profile of nucleoside reverse transcriptase inhibitor (NRTI) BMS-986001 against known NRTI resistance mutations. *Antimicrob. Agents Chemother.* **57(11)**, 5500-5508 (2013).

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, 10/04/2023

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