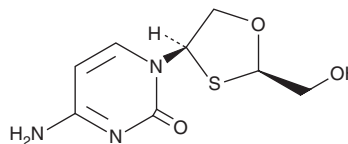


PRODUCT INFORMATION



Apricitabine Item No. 36037

CAS Registry No.: 160707-69-7
Formal Name: 4-amino-1-[(2R,4R)-2-(hydroxymethyl)-1,3-oxathiolan-4-yl]-2(1H)-pyrimidinone
Synonyms: AVX754, BCH 10618, (-)-BCH 10652, (-)-dOTC, SPD-754
MF: C₈H₁₁N₃O₃S
FW: 229.3
Purity: ≥95%
UV/Vis.: λ_{max}: 277 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

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

Description

Apricitabine is a nucleoside reverse transcriptase inhibitor (NRTI) and derivative of 2'-deoxycytidine (Item Nos. 34708 | 30125).¹ It is phosphorylated intracellularly to its active form, apricitabine-5'-triphosphate, which competes with a monophosphate form for binding to HIV-1 reverse transcriptase and incorporation into the nascent viral DNA, inducing chain termination and HIV-1 replication. Apricitabine reduces HIV-1 infectivity in MT-4 and U937 cells (IC₅₀s = 2.8 and 0.4 μM, respectively), as well as in isolated human cord blood mononuclear cells (CBMCs; IC₅₀ = 0.4 μM).²

References

1. Cox, S. and Southby, J. Apricitabine--a novel nucleoside reverse transcriptase inhibitor for the treatment of HIV infection that is refractory to existing drugs. *Expert Opin. Investig. Drugs* **18(2)**, 199-209 (2009).
2. Mansour, T.S., Jin, H., Wang, W., et al. Structure-activity relationships among a new class of antiviral heterosubstituted 2',3'-dideoxynucleoside analogues. *Nucleos. Nucleot.* **14(3-5)**, 627-635 (1995).

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, 12/19/2022

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