

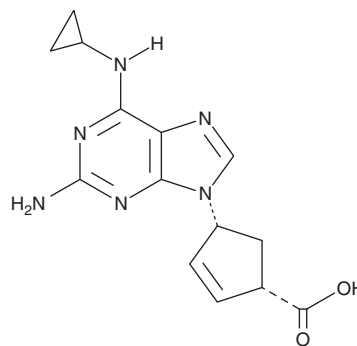
PRODUCT INFORMATION



Abacavir Carboxylate

Item No. 27777

CAS Registry No.: 384380-52-3
Formal Name: (1S,4R)-4-[2-amino-6-(cyclopropylamino)-9H-purin-9-yl]-2-cyclopentene-1-carboxylic acid
MF: C₁₄H₁₆N₆O₂
FW: 300.3
Purity: ≥95%
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

Abacavir carboxylate is supplied as a solid. A stock solution may be made by dissolving the abacavir carboxylate in the solvent of choice, which should be purged with an inert gas. Abacavir carboxylate is soluble in organic solvents such as DMSO and methanol.

Description

Abacavir carboxylate is an inactive metabolite of the HIV-1 reverse transcriptase inhibitor abacavir (Item No. 14746).^{1,2} It is formed from abacavir *via* reactive aldehyde intermediates that can form adducts with proteins on valine residues.²

References

1. Martin, M.A. and Kroetz, D.L. Abacavir pharmacogenetics--from initial reports to standard of care. *Pharmacotherapy* **33(7)**, 765-775 (2013).
2. Charneira, C., Godinho, A.L.A., Oliveira, M.C., *et al.* Reactive aldehyde metabolites from the anti-HIV drug abacavir: Amino acid adducts as possible factors in abacavir toxicity. *Chem. Res. Toxicol.* **24(12)**, 2129-21241 (2011).

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, 12/22/2022

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