

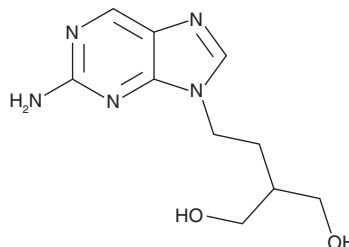
PRODUCT INFORMATION



6-Deoxypenciclovir

Item No. 35602

CAS Registry No.: 104227-86-3
Formal Name: 2-[2-(2-amino-9H-purin-9-yl)ethyl]-1,3-propanediol
Synonym: BRL 42359
MF: C₁₀H₁₅N₅O₂
FW: 237.3
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

6-Deoxypenciclovir is supplied as a solid. A stock solution may be made by dissolving the 6-deoxypenciclovir in the solvent of choice, which should be purged with an inert gas. 6-Deoxypenciclovir is slightly soluble in methanol and DMSO.

Description

6-Deoxypenciclovir is an inactive metabolite of the antiviral prodrug famciclovir (Item No. 23834).^{1,2} It is formed from famciclovir via hydrolysis and is further metabolized via oxidation by aldehyde oxidase to the active metabolite penciclovir (Item No. 22918). 6-Deoxypenciclovir has been used in the synthesis of prodrug forms of penciclovir.³

References

1. Rashidi, M.R., Smith, J.A., Clarke, S.E., *et al.* In vitro oxidation of famciclovir and 6-deoxypenciclovir by aldehyde oxidase from human, guinea pig, rabbit, and rat liver. *Drug Metab. Dispos.* **25**(7), 805-813 (1997).
2. Groth, A.D., Contreras, M.T., Kado-Fong, H.K., *et al.* In vitro cytotoxicity and antiviral efficacy against feline herpesvirus type 1 of famciclovir and its metabolites. *Vet. Ophthalmol.* **17**(4), 268-274 (2014).
3. Kim, D.K., Lee, N., Ryu, D.H., *et al.* Synthesis and evaluation of 2-amino-9-(3-acyloxymethyl-4-alkoxycarbonyloxybut-1-yl)purines and 2-amino-9-(3-alkoxycarbonyloxymethyl-4-alkoxycarbonyloxybut-1-yl)purines as potential prodrugs of penciclovir. *Bioorg. Med. Chem.* **7**(8), 1715-1725 (1999).

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, 12/12/2022

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