

PRODUCT INFORMATION



PC-766B

Item No. 20587

CAS Registry No.: 108375-77-5
Formal Name: 16-[(1S,2R,3S)-3-[(2R,4R,5S,6R)-4-[(2,6-dideoxy-β-D-arabino-hexopyranosyl)oxy] tetrahydro-2-hydroxy-5-methyl-6-[(1E,3E)-1,3-pentadien-1-yl]-2H-pyran-2-yl]-2-hydroxy-1-methylbutyl]-8-hydroxy-3,15-dimethoxy-5,7,9,11-tetramethyl-(3Z,5E,7R,8S,9S,11E,13E,15S,16R)-oxacyclohexadeca-3,5,11,13-tetraen-2-one

Synonym: Antibiotic PC-766B

MF: C₄₃H₆₈O₁₂

FW: 777.0

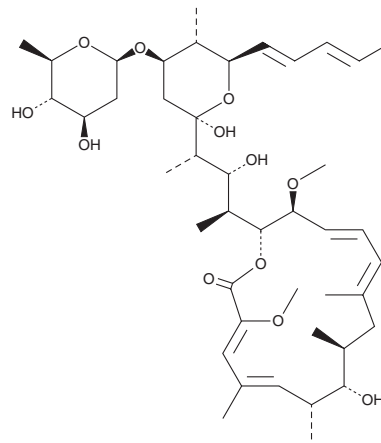
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Nocardia brasiliensis*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PC-766B is supplied as a solid. A stock solution may be made by dissolving the PC-766B in the solvent of choice, which should be purged with an inert gas. PC-766B is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

PC-766B is a macrolide antibiotic produced by *N. brasiliensis* that is active against Gram-positive bacteria and some fungi, yet inactive against Gram-negative bacteria.¹ It also demonstrates antitumor activity against mouse P388 leukemia and B16 melanoma cells *in vitro* (IC₅₀s = 0.1 and 0.4 ng/ml, respectively) and *in vivo*, but weakly inhibits the Na⁺,K⁺-ATPase compared to structurally related antibiotics.¹

Reference

1. Kumagai, K., Taya, K., Fukui, A., *et al.* PC-766B, a new macrolide antibiotic produced by *Nocardia brasiliensis*. I. Taxonomy, fermentation and biological activity. *J. Antibiot. (Tokyo)* **46(6)**, 972-978 (1993).

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, 10/18/2022

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