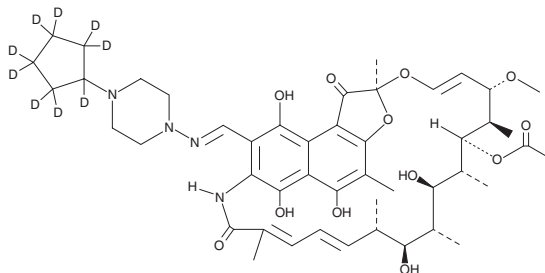


PRODUCT INFORMATION



Rifapentine-d₉ Item No. 28708

Formal Name: 3-[[[(4-cyclopentyl-1-piperazinyl)imino]methyl]-rifamycin-d₉
Synonym: Cyclopentylrifampicin-d₉
MF: C₄₇H₅₅D₉N₄O₁₂
FW: 886.1
Chemical Purity: ≥98% (Rifapentine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
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

Rifapentine-d₉ is intended for use as an internal standard for the quantification of rifapentine (Item No. 20307) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Rifapentine-d₉ is supplied as a solid. A stock solution may be made by dissolving the rifapentine-d₉ in the solvent of choice, which should be purged with an inert gas. Rifapentine-d₉ is soluble in the organic solvents DMSO, methanol, and acetonitrile.

Description

Rifapentine is a long-lasting rifamycin antibiotic that inhibits DNA-dependent RNA polymerases and is effective against *Mycobacterium* species, including *M. tuberculosis*, as well as experimental infections caused by *S. aureus*, *S. pyogenes* group A, *S. pneumoniae*, and *K. pneumoniae*.^{1,2}

References

1. Arioli, V., Berti, M., Carniti, G., *et al.* Antibacterial activity of DL 473, a new semisynthetic rifamycin derivative. *J. Antibiot. (Tokyo)* **34(8)**, 1026-1032 (1981).
2. Sharma, S.K., Sharma, A., Kadiravan, T., *et al.* Rifamycins (rifampicin, rifabutin and rifapentine) compared to isoniazid for preventing tuberculosis in HIV-negative people at risk of active TB. *Cochrane Database Syst. Rev.* **7**, CD007545 (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, 10/13/2022

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