

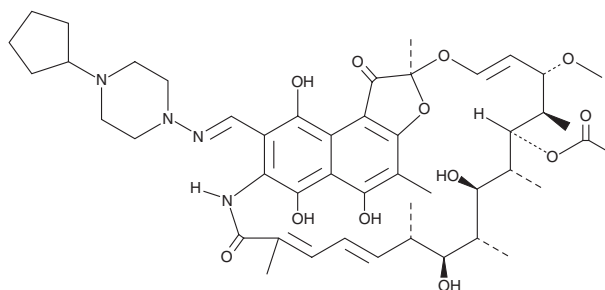
# PRODUCT INFORMATION



## Rifapentine

Item No. 20307

**CAS Registry No.:** 61379-65-5  
**Formal Name:** 3-[[[4-(cyclopentyl-1-piperazinyl) imino]methyl]-rifamycin  
**Synonyms:** Cyclopentylrifampicin, DL-473, MDL 473,  
**MF:** C<sub>47</sub>H<sub>64</sub>N<sub>4</sub>O<sub>12</sub>  
**FW:** 877.0  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 237, 253, 336, 478 nm  
**Supplied as:** A crystalline solid  
**Storage:** Room temperature  
**Stability:** ≥4 years



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

### Laboratory Procedures

Rifapentine is supplied as a crystalline solid. A stock solution may be made by dissolving the rifapentine in the solvent of choice, which should be purged with an inert gas. Rifapentine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of rifapentine in ethanol is approximately 2 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Rifapentine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, rifapentine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Rifapentine has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### 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*.<sup>1,2</sup>

### 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**, (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, 12/01/2022

#### 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