

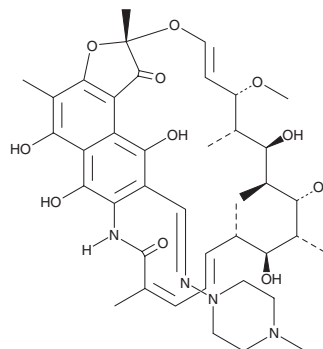
PRODUCT INFORMATION



25-Desacetyl Rifampicin

Item No. 34817

CAS Registry No.: 16783-99-6
Formal Name: 25-O-deacetyl-3-[(E)-[4-methyl-1-piperazinyl]imino]methyl]-rifamycin
Synonym: DesacetylRifampicin
MF: C₄₁H₅₆N₄O₁₁
FW: 780.9
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

25-Desacetyl rifampicin is supplied as a solid. A stock solution may be made by dissolving the 25-Desacetyl rifampicin in the solvent of choice, which should be purged with an inert gas. 25-Desacetyl rifampicin is slightly soluble in DMSO.

Description

25-Desacetyl rifampicin is a major active metabolite of the rifamycin antibiotic rifampicin (Item No. 14423).¹ 25-Desacetyl rifampicin is active against *M. smegmatis* (MIC₉₉ = 2.66 μM).

Reference

1. Kigonde, E.M., Njoroge, M., Singh, K., *et al.* Synthesis and synergistic antimycobacterial screening of chlorpromazine and its metabolites. *Med. Chem. Commun.* 5, 502-506 (2014).

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, 12/13/2022

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