

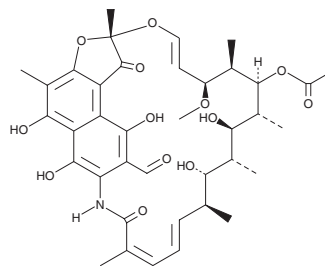
PRODUCT INFORMATION



3-formyl Rifamycin

Item No. 20260

CAS Registry No.: 13292-22-3
Formal Name: 3-formyl-rifamycin
Synonyms: 3-Formyl Rifamycin SV,
NCI 145-635, Rifaldehyde
MF: C₃₈H₄₇NO₁₃
FW: 725.8
Purity: ≥95%
UV/Vis.: λ_{max}: 220, 240, 262, 324, 489 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

3-formyl Rifamycin is an intermediate of rifampicin (Item No. 14423) that has been used to develop several rifamycin derivatives with antibiotic activity that targets RNA synthesis.¹

Reference

1. Riva, S. and L. G. Silvestri. Rifamycins: A general view. *Annu. Rev. Microbiol.* **26**, 199-224 (1982).

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, 11/30/2022

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