

Rifampicin

Bacterial DNA-dependent RNA polymerase inhibitor

A selective inhibitor of bacterial DNA-dependent RNA polymerase (RNAP). Has no effect on mammalian RNA polymerase. The RNAP-rifampicin complex locks the promoter in an abortive initiation reaction making short oligoribonucleotides which diffuse out of the active site. Inhibits α -synuclein fibrillation and disaggregates fibrils. Potent inducer of hepatic cytochrome P450 enzymes.

Citations: 9

[View Online »](#)

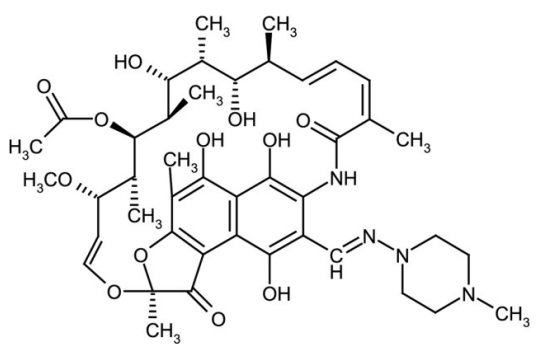
Ordering Information

[Order Online »](#)

ALX-380-071-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Rifampin, 3-(4-Methylpiperazinyl-iminomethyl)rifamycin SV

Appearance Orange to brown powder.

CAS 13292-46-1

Couple Target Bacterial DNA-dependent RNA polymerase

Couple Type Inhibitor

Formula $C_{43}H_{58}N_4O_{12}$

MI 14: 8216

MW 823.0

Purity ≥97% (Assay)

RTECS VJ7000000

Solubility Soluble in DMSO (25mg/ml) or methanol (25mg/ml).

Source Semisynthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-](mailto:info-usa@enzolifesciences.com)
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-](mailto:info-be@enzolifesciences.com)
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-](mailto:info-fr@enzolifesciences.com)
be@enzolifesciences.com

France
Phone: +33 472 440 655
[info-](mailto:info-fr@enzolifesciences.com)
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
[info-](mailto:info-de@enzolifesciences.com)
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-](mailto:info-uk@enzolifesciences.com)
uk@enzolifesciences.com