

Enniatin B

Ionophore antibiotic

Ionophore antibiotic. Inhibitor of the *S. cerevisiae* ABC transporter Pdr5p. Inhibitor of acyl-CoA-cholesterol acyltransferase and of phosphodiesterase. Potential anticancer compound.

Citations: 9

[View Online »](#)

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

Appearance White to off-white crystalline solid.

CAS 917-13-5

Couple Target ACAT, MDR, PDE

Couple Type Inhibitor

Formula $C_{33}H_{57}N_3O_9$

MI **14: 3585**

MW 639.8

Purity ≥90% (HPLC)

Solubility Soluble in DMSO or other organic solvents; almost insoluble in water.

Source Isolated from *Fusarium orthoceras var. enniatum*.

Last modified: May 29, 2024



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

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

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

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

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

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