

Ascomycin

Immunosuppressant

Ascomycin is a close structural analog of FK506 and is a potent immunosuppressant. It binds to FKBP (FK506-binding protein) and suppresses its proline rotamase activity. The ascomycin-FKBP complex inhibits calcineurin, a type 2B phosphatase.

Citations: 8

[View Online »](#)

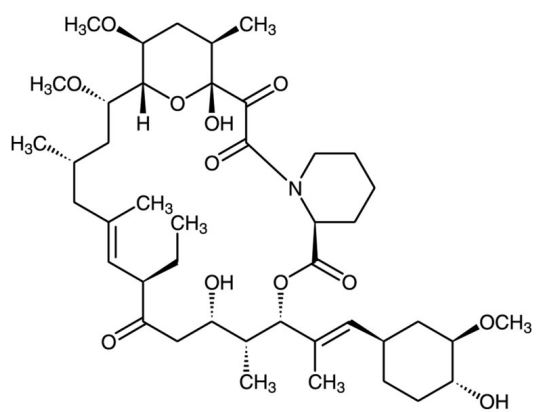
Ordering Information

[Order Online »](#)

BML-A165-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Immunomycin
Appearance	Off-white solid
CAS	104987-12-4; also registered as 11011-38-4
Couple Target	FKBP, Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	$C_{43}H_{69}NO_{12}$
MW	792
Purity	≥90% (HPLC, TLC)
Solubility	Soluble in DMSO, methanol or ethanol.
Source	Isolated from <i>Streptomyces</i> sp.



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-
fr@enzolifesciences.com](mailto:info-fr@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)