

Tunicamycin

GlcNAc phosphotransferase inhibitor

Nucleoside antibiotic. Inhibits protein N-glycosylation by blocking the transfer of N-acetylglucosamine 1-phosphate to dolichol monophosphate. Induces ER-stress following the inhibition of N-linked glycosylation. Arrests cell cycle in late G1. Unstable at acidic pH.

Citations: 28

[View Online »](#)

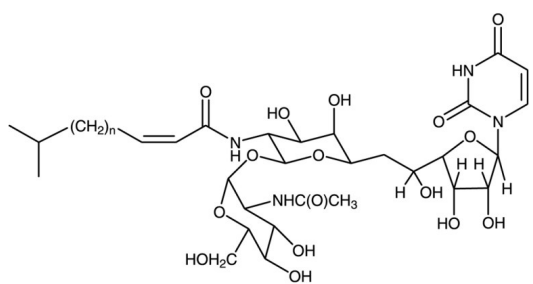
Ordering Information

[Order Online »](#)

BML-CC104-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid to tan solid.
CAS	11089-65-9
Couple Target	GlcNAc phosphotransferase
Couple Type	Inhibitor
Formula	$C_{38}H_{62}N_4O_{16}$ (Tun B)
MI	14: 9819
MW	Avg. 840.0
Purity	≥99% (TLC)
Purity Detail	Mixture of tunicamycin A, B, C and D.
RTECS	YO7980200
Solubility	Soluble in DMSO (>10mg/ml) or methanol (5mg/ml; warm).
Source	Isolated from <i>Streptomyces lysosuperficus</i> .

Technical Info / Product Notes



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Last modified: May 29, 2024

Replacement for **ADI-908-297**.