

Thiostrepton

Elongation factor G inhibitor

Thiazole-containing peptide antibiotic. Inhibits protein synthesis by preventing binding of GTP to the 50S ribosomal subunit. Inhibits binding of elongation factor G (EF-G) and the dissociation of EF-G from the ribosome.

Citations: 12

[View Online »](#)

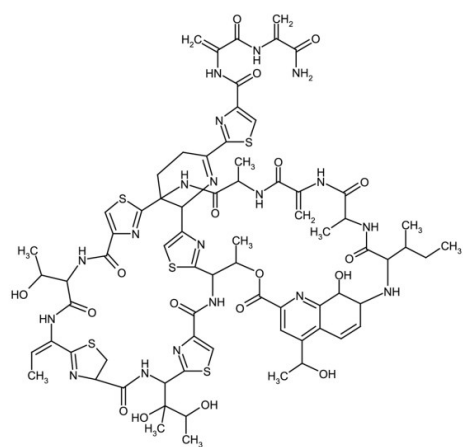
Ordering Information

[Order Online »](#)

ALX-380-261-G001	1g
ALX-380-261-G005	5g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Activity	≥900 U/mg
Appearance	White to off-white solid.
CAS	1393-48-2
Couple Target	Elongation factor, Ribosome
Couple Type	Inhibitor, Ligand
Formula	$C_{72}H_{85}N_{19}O_{18}S_5$
MI	14: 9364
MW	1664.9
RTECS	XN6300100
Solubility	Soluble in acetic acid, DMSO (10mg/ml) or dimethyl formamide; insoluble in water.
Source	Isolated from <i>Streptomyces azureus</i> .

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-
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)