

Fusidic acid . sodium salt

Elongation factor G inhibitor

Antibiotic which suppresses nitric oxide (NO) toxicity in pancreatic islet cells. Inhibits protein synthesis by inhibition of elongation factor G at the level of the ribosome.

Citations: 7

[View Online »](#)

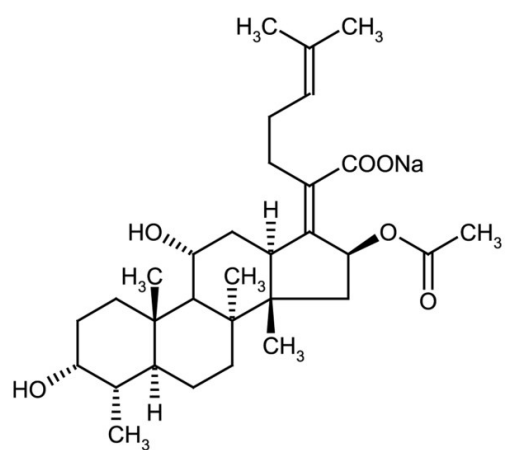
Ordering Information

[Order Online »](#)

ALX-380-011-M050	50mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white solid.

CAS 751-94-0

Couple Target Elongation factor

Couple Type Inhibitor

Formula $C_{31}H_{47}O_6 \cdot Na$

MI **14:** 4317

MW 515.7 . 23.0

Purity ≥98% (TLC)

Solubility Soluble in water (50mg/ml).

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)