

Reveromycin A

Cell cycle inhibitor

Antibiotic. Dominant analog of the reveromycin complex. Inhibitor of the mitogenic activity of epidermal growth factor (EGF). G1 phase cell cycle inhibitor, selectively inhibiting isoleucyl-tRNA synthetase. Displays antiproliferative as well as potent antifungal activity. Induces apoptosis in osteoclasts thus inhibiting bone resorption.

Citations: 8

[View Online »](#)

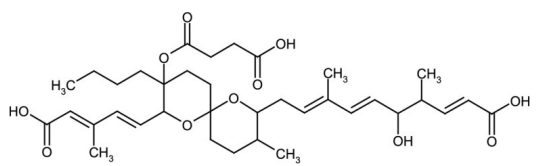
Ordering Information

[Order Online »](#)

ALX-380-216-MC01	0.1mg
ALX-380-216-MC25	0.25mg

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 tan solid.

CAS 134615-37-5

Couple Target EGF, Isoleucyl-tRNA synthetase

Couple Type Inhibitor

Formula $C_{36}H_{52}O_{11}$

Formulation Lyophilized.

MW 660.8

Purity ≥99% (HPLC)

Solubility Soluble in ethyl acetate, 100% ethanol, methanol, dimethyl formamide or DMSO.

Source Isolated from *Streptomyces* sp. MST-AS5344.

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)