

Fumitremorgin C

BCRP inhibitor

Mycotoxin. Tremorgenic. Potent and specific inhibitor of the breast cancer resistance protein (BCRP; ABCG2). Reverses multidrug resistance mediated by BCRP and increases cytotoxicity of several anticancer agents *in vitro*.

Citations: 17

[View Online »](#)

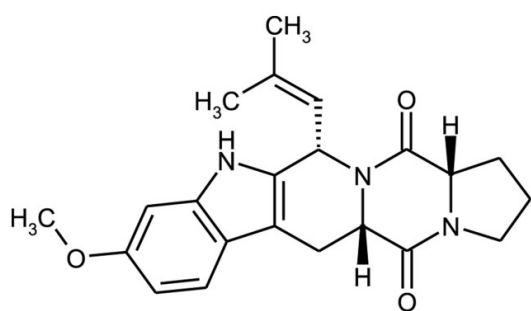
Ordering Information

[Order Online »](#)

ALX-350-127-C250	250µg
------------------	-------

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

Alternative Name FTC, SM-Q

Appearance White to light brown powder.

CAS 118974-02-0

Couple Target BCRP

Couple Type Inhibitor

Formula $C_{22}H_{25}N_3O_3$

Identity Determined by ¹H-NMR.

MW 379.5

Purity ≥96% (HPLC)

Solubility Soluble in DMSO, methanol or chloroform.

Source Isolated from *Aspergillus fumigatus*.

Technical Info / Product Notes Optical Rotation: $[\alpha]_D = +19^\circ \pm 1^\circ$ (c=1.0 in methanol).

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

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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