

Tryprostatin A

Cell cycle blocker and BCRP inhibitor

Diketopiperazine. Arrests cell cycle progression at the G2/M phase. Inhibitor of MAP-dependent microtubule assembly. Inhibits the ABC-transporter breast cancer resistance protein (BCRP) and thereby reverses the resistance of cancer cells against antitumor drugs.

Citations: 5

[View Online »](#)

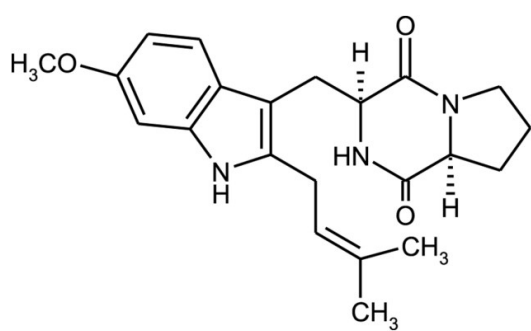
Ordering Information

[Order Online »](#)

ALX-380-090-C500	500µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White powder.
CAS	171864-80-5
Couple Target	BCRP
Couple Type	Inhibitor
Formula	$C_{22}H_{27}N_3O_3$
MW	381.5
Purity	≥90% (HPLC)
Solubility	Soluble in methanol, 100% ethanol or DMSO.
Source	Isolated from <i>Aspergillus fumigatus</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)