

Atpenin A5 (synthetic)

Mitochondrial complex II inhibitor

Antibiotic. Potent and specific inhibitor of mitochondrial complex II (succinate-ubiquinone oxidoreductase).

Citations: 13

[View Online »](#)

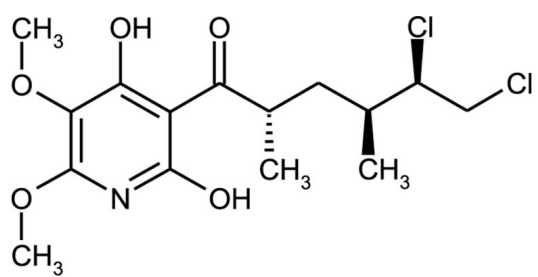
Ordering Information

[Order Online »](#)

ALX-380-313-MC25	0.25mg
ALX-380-313-M001	1mg

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

Alternative Name	3-[(2S,4S,5R)-5,6-Dichloro-2,4-dimethyl-1-oxohexyl]-4-hydroxy-5,6-dimethoxy-2(1H)-pyridinone
Appearance	White to off-white powder.
CAS	119509-24-9
Couple Target	Mitochondrial complex
Couple Type	Inhibitor
Formula	C ₁₅ H ₂₁ Cl ₂ NO ₅
MW	366.2
Purity	≥95% (HPLC)
RTECS	CJ8800000
Solubility	Soluble in acetone, acetonitrile, chloroform, ethyl acetate, DMSO, methanol or 100% ethanol; insoluble in water or hexane.
Source	Synthetic. Originally isolated from <i>Penicillium</i> sp. strain FO-125.

The IC₅₀ value against bovine heart complex II is 3.6nM (which is ~300-fold lower than the IC₅₀ value of carboxin (1.1μM)). It also inhibits fumarate reductase of *Ascaris suum* (IC₅₀ = 12nM). Inhibition of *E. coli* succinate dehydrogenase is less potent (IC₅₀ = 5μM). By co-crystallization studies of atpenin A5 and complex II, the binding site of atpenin A5 was found to be the quinone-binding site of complex II. Additionally, atpenin A5 has been shown to have a protective action against ischemia-reperfusion via the activation of mitochondrial K_{ATP} channels.

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)