

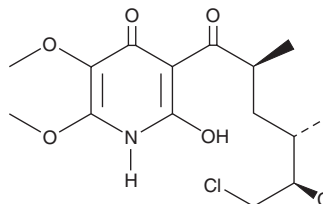
PRODUCT INFORMATION



Atpenin A5

Item No. 11898

CAS Registry No.: 119509-24-9
Formal Name: 3-[(2S,4S,5R)-5,6-dichloro-2,4-dimethyl-1-oxohexyl]-4-hydroxy-5,6-dimethoxy-2(1H)-pyridinone
Synonym: Antibiotic FO-125A5
MF: C₁₅H₂₁Cl₂NO₅
FW: 366.2
Purity: ≥95%
UV/Vis.: λ_{max}: 237, 272, 324 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic; originally from fungal strain FO-125



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Atpenin A5 is supplied as a crystalline solid. A stock solution may be made by dissolving the atpenin A5 in the solvent of choice. Atpenin A5 is soluble in organic solvents such as ethanol, methanol, and DMSO, which should be purged with an inert gas.

Description

Mitochondrial complex II (succinate dehydrogenase or succinate:ubiquinone oxidoreductase) is a functional member of the Krebs cycle and the aerobic respiratory chain that couples the oxidation of succinate to fumarate with the reduction of quinone to quinol. Atpenin A5, an antifungal antibiotic isolated from *Penicillium* sp. found in soil, is a highly specific ubiquinone-binding site inhibitor of succinate dehydrogenase (IC₅₀s = 12 and 3.7 nM in nematode and mammalian mitochondria, respectively, versus IC₅₀s > 100 μM for inhibition of complex I and complex III enzymes).^{1,2} Atpenin A5 has cardioprotective effects against simulated ischemia-reperfusion injury in cardiomyocytes. Several mechanisms through which this occurs, including activation of mitochondrial ATP-sensitive potassium channels or modulation of mitochondrial reactive oxygen species generation, have been proposed.^{3,4}

References

1. Horsefield, R., Yankovskaya, V., Sexton, G., *et al.* Structural and computational analysis of the quinone-binding site of complex II (succinate-ubiquinone oxidoreductase): A mechanism of electron transfer and proton conduction during ubiquinone reduction. *J. Biol. Chem.* **281**(11), 7309-7316 (2006).
2. Miyadera, H., Shiomi, K., Ui, H., *et al.* Atpenins, potent and specific inhibitors of mitochondrial complex II (succinate-ubiquinone oxidoreductase). *Proc. Natl. Acad. Sci. USA* **100**(2), 473-477 (2003).
3. Wojtovich, A.P. and Brookes, P.S. The complex II inhibitor atpenin A5 protects against cardiac ischemia-reperfusion injury via activation of mitochondrial KATP channels. *Basic Res. Cardiol.* **104**(2), 121-129 (2009).
4. Dröse, S., Bleier, L., and Brandt, U. A common mechanism links differently acting complex II inhibitors to cardioprotection: Modulation of mitochondrial reactive oxygen species production. *Mol. Pharmacol.* **79**(5), 814-822 (2011).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/13/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM