

Sanguinarine chloride

Na⁺/K⁺- and Mg²⁺-ATPase inhibitor

Na⁺/K⁺-and Mg²⁺-ATPase inhibitor. Antimicrobial, anti-inflammatory and antioxidant properties. Induces apoptosis. Modulates the Bcl-2 family. Shows antiplaque activity in humans.

Citations: 9

[View Online »](#)

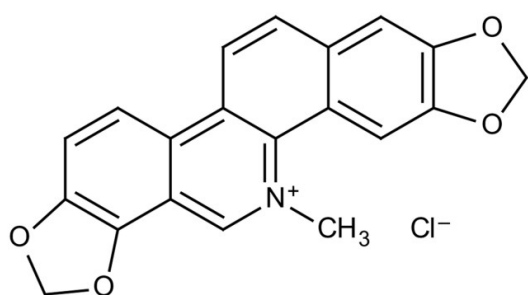
Ordering Information

[Order Online »](#)

ALX-350-076-M005	5mg
------------------	-----

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 Pseudochelerythrine

Appearance Orange solid.

CAS 5578-73-4

Couple Target ATPase

Couple Type Inhibitor

Formula $C_{20}H_{14}NO_4Cl$

MI 14: 8356

MW 367.8

Purity ≥98% (HPLC)

RTECS VP5220000

Solubility Soluble in methanol; slightly soluble in water.

Source Isolated from the leaves and stems of *Macleaya cordata* and *M. microcarpa*.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)