

Resibufogenin

Na⁺/K⁺ ATPase inhibitor

Specific Na⁺/K⁺-ATPase inhibitor.

Citations: 2

[View Online »](#)

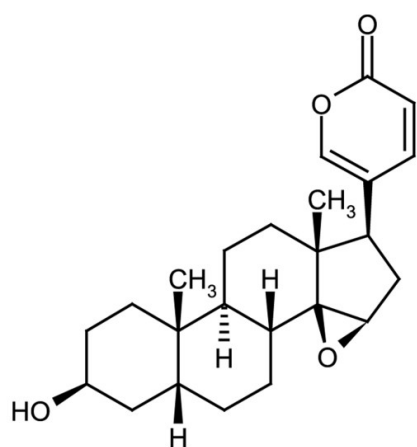
Ordering Information

[Order Online »](#)

ALX-350-286-M010	10mg
------------------	------

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 (3β,5β,15β)-14,15-Epoxy-3-hydroxy-5-bufa-20,22-dienolide

Appearance White to off-white solid.

CAS 465-39-4

Couple Target ATPase

Couple Type Inhibitor

Formula $C_{24}H_{32}O_4$

Identity Determined by MS.

MI **14:** 8146

MW 384.5

Purity ≥98% (HPLC)

RTECS EI2975000

Solubility Soluble in chloroform.

Source Isolated from *Bufo* venom.



ENZO LIFE SCIENCES,
INC.
May 29, 2024
info-
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
be@enzolifesciences.com](mailto:info-be@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)