

Ouabain . octahydrate

Na⁺/K⁺ ATPase inhibitor

A steroidal inhibitor of cardiac sodium pumps (Na⁺/K⁺ ATPase) which has been recognized as an endogenous adrenal hormone. In the CNS it can be used as a tool to induce in vivo excitotoxicity. Ouabain, as well as other endogenous digitalis-like compounds, can down-regulate the expression of 14-3-3 proteins in rat lens.

Citations: 16

[View Online »](#)

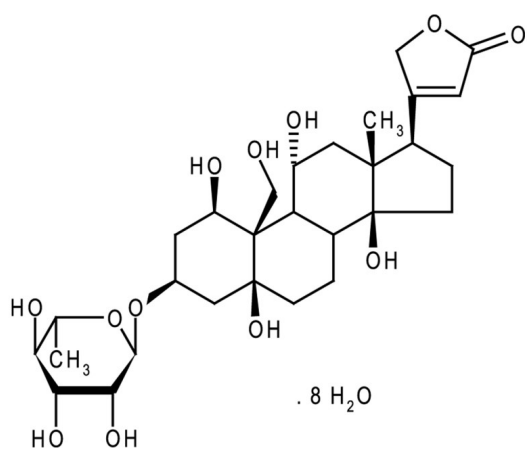
Ordering Information

[Order Online »](#)

ALX-350-066-M100	100mg
------------------	-------

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	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Strophanthin G, 3-[(6-Deoxy- α -L-mannopyranosyl)oxy]-1,5,11 α -14,19-penta-hydroxycard-20(22)-enolide
Appearance	White solid.
CAS	11018-89-6
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{29}H_{44}O_{12} \cdot 8H_2O$
MI	14: 6901
MW	584.6 . 144.2
Purity	$\geq 98\%$ (TLC)
Solubility	Soluble in water (>25mg/ml).

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)