

Ryanodine (high purity)

Ca²⁺ channel activator

Induces the release of intracellular Ca²⁺. Ryanodine receptors mediate adenine nucleotide, Ca²⁺ and caffeine-sensitive release of Ca²⁺ from sarcoplasmic reticulum of cardiac and skeletal muscle and may also mediate intracellular Ca²⁺ release in neurons. Biphasic effect on sarcoplasmic reticulum calcium channels; Nanomolar levels open the channel and micromolar levels close it.

Citations: 27

[View Online »](#)

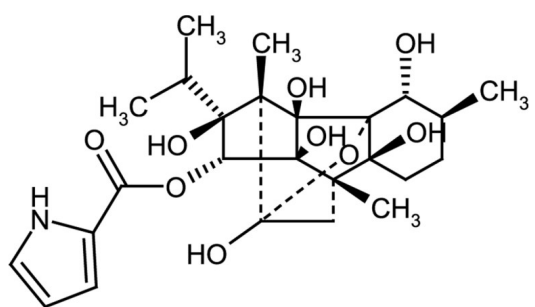
Ordering Information

[Order Online »](#)

ALX-630-062-M001	1mg
ALX-630-062-M005	5mg

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 -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white solid.

CAS 15662-33-6

Couple Target Calcium channel

Couple Type Activator

Formula $C_{25}H_{35}NO_9$

MI 14: 8307

MW 493.5

Purity ≥99% (HPLC)

RTECS VM4025000

Solubility Soluble in methanol or DMSO (>29mg/ml), in 100% ethanol (16mg/ml); slightly soluble in water (warm).

Source Isolated from *Ryania speciosa*, *Flacourtiaceae*.



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)