

Hellebrin

Na⁺/K⁺ ATPase inhibitor

Cardiotonic glycoside which presents a bufadienolide-steroid structure. Besides ouabain (Prod. No. ALX-350-066), from the class of cardenolides hellebrin is the second unique water soluble cardiotonic glycoside. Extremely potent inhibitor of Na⁺/K⁺-ATPase (sodium pump) blocking the active efflux of Na⁺ and reuptake of K⁺ of this membrane enzyme. Shows a positive inotropic effect by increasing the intracellular Ca²⁺ concentration. Strong immunosuppressor inhibiting T cell activity with much higher potency than cortisol or cyclosporin A (Prod. No. BML-A195). Induces caspase-dependent apoptosis.

Citations: 5

[View Online »](#)

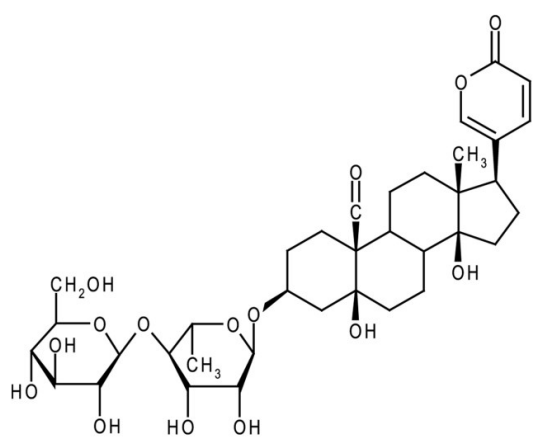
Ordering Information

[Order Online »](#)

ALX-350-105-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 -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 3β-(O4-β-D-Glucopyranosyl-α-L-rhamnopyranosyloxy)-5,14-dihydroxy-19-oxo-5β,14-β-bufa-20,22-dienolide

Appearance White to off-white solid.

CAS 13289-18-4

Couple Target ATPase

Couple Type Inhibitor

Formula C₃₆H₅₂O₁₅

MW 724.8

Purity ≥99% (HPLC)

Solubility Soluble in water (4.4mg/ml).

Source Isolated from *Helleborus sp.*

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)