

Bufotalin

Na⁺/K⁺ ATPase inhibitor

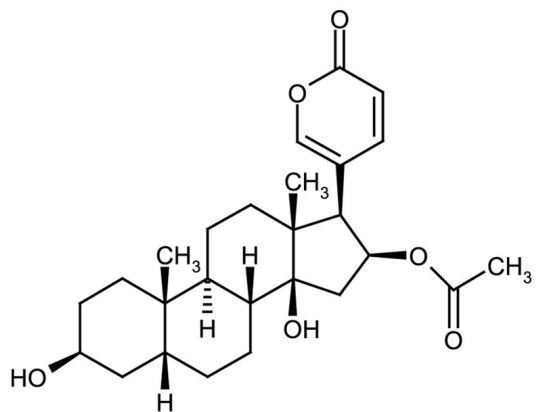
Less potent Na⁺/K⁺-ATPase inhibitor than other bufadienolides. Potent immunosuppressor.

Citations: 2

[View Online »](#)

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (3β,5β,16β)-16-(Acetyloxy)-3-,14-dihydroxybufa-20,22-dienolide, 3β,14,16β-Trihydroxy-5β bufa-20,22-dienolide 16-acetate

Appearance White to off-white solid.

CAS 471-95-4

Couple Target ATPase

Couple Type Inhibitor

Formula $C_{26}H_{36}O_6$

MI **14:** 1477

MW 444.6

Purity ≥98% (HPLC)

Solubility Soluble in 100% ethanol or chloroform.

Source Isolated from *Bufois* venom.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com