

Bufalin

ATPase inhibitor

Cardiotonic steroid specific Na^+/K^+ -ATPase inhibitor. More potent than ouabain (Prod. No. BML-CM109). Induces cell differentiation. Induces apoptosis in U937 and HL-60 cells by activation of AP-1 via a MAPK pathway that involves JNK. Regulates Tiam1 gene. Down regulates expression of 14-3-3 protein.

Citations: 11

[View Online »](#)

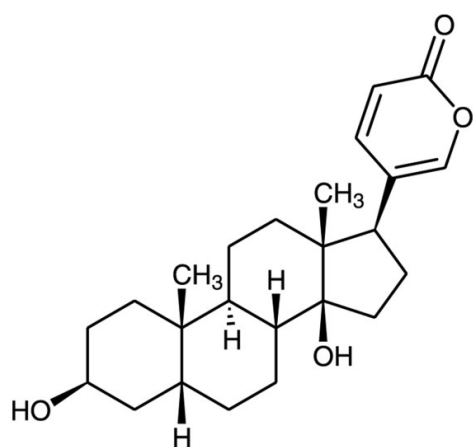
Ordering Information

[Order Online »](#)

BML-AP303-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(3β,5β)-3,14-Dihydroxybufa-20,22-dienolide, 5β,20(22)-Bufadienolide-3β,14-diol
Appearance	White to off-white solid.
CAS	465-21-4
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{24}H_{34}O_4$
MI	14: 1471
MW	386.5
Purity	≥97% (HPLC)
RTECS	EI2962500
Solubility	Soluble in DMSO (25 mg/ml) or 100% ethanol (25 mg/ml, warm).

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-fr@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