

# Novobiocin . sodium salt

## HSP90 inhibitor

Antibiotic. Inhibitor of HSP90. Interacts with the C-terminal ATP-binding domain of HSP90 in contrast to the benzoquinone ansamycins 17-AAG (Prod. No. BML-EI308), geldanamycin (Prod. No. BML-EI280), herbimycin A (Prod. No. BML-EI227) and the chemically unrelated radicicol (Prod. No. BML-EI285), which bind to the N-terminal ATP-binding site of HSP90. Inhibitor of DNA gyrase. Useful for the production of positively supercoiled plasma DNA, targeting the nucleotide-binding site of gyrase B. Inhibits retrovirus RNA-dependent DNA polymerase. Potent inhibitor of ADP ribosylation. Inhibits LPS-induced production of pro-inflammatory cytokines, such as TNF- $\alpha$ , IL-1, IL-6 and IL-10. Inhibits protein synthesis and alters the phosphorylation state of several cytosolic proteins. Reverses etoposide resistance in non-P-glycoprotein expressing multidrug resistant (MDR) tumor cell lines. Inducer of CD38 on cells of the myelomonocytic lineage. Forms ion channels in lipid bilayers. Inhibitor of breast cancer resistance protein (BCRP)

Citations: 29

[View Online »](#)

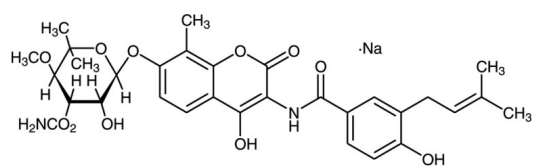
## Ordering Information

[Order Online »](#)

BML-A256-0001	1g
BML-A256-0005	5g

## 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

Activity	~900 µg/mg
Alternative Name	Albamycin, Streptonivicin, N-[7-[[3-O-Aminocarbonyl-6-deoxy-5-C-methyl-4-O-β-L-lyxo-hexapyranosyl]oxy]-4-hydroxy-8-methyl-2-oxo-2H-1-benzopyran-3-yl]-4-hydroxy-3-(3-methyl-2-butenyl)benzamide
Appearance	White to off-white powder.
CAS	1476-53-5
Couple Target	HSP90
Couple Type	Inhibitor
Excitation Maximum	307nm
Formula	$C_{31}H_{35}N_2O_{11} \cdot Na$
MI	14: 6722
MW	634.6
RTECS	RD5425000

## Solubility

Soluble in water (100mg/ml), ethanol or DMSO.



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)

Last modified: May 29, 2024