

# Valinomycin

## Potassium ionophore

A potassium ionophore. Induces mitochondrial swelling. Induces apoptosis in CHO cells and other cell lines. Antagonizes endothelin-induced vasoconstriction ( $IC_{50}=0.3\mu M$ ).

**Citations:** 18

[View Online »](#)

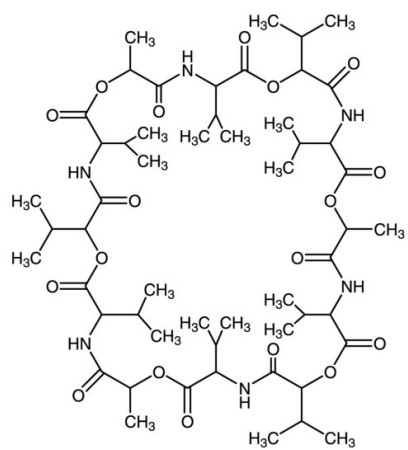
## Ordering Information

[Order Online »](#)

|                |      |
|----------------|------|
| BML-KC140-0025 | 25mg |
|----------------|------|

## Manuals, SDS & CoFA

[View Online »](#)



## Handling & Storage

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

## Regulatory Status

RUO - Research Use Only

## Product Details

|            |                                                                            |
|------------|----------------------------------------------------------------------------|
| Appearance | White crystalline solid.                                                   |
| CAS        | 2001-95-8                                                                  |
| Formula    | $C_{54}H_{90}N_6O_{18}$                                                    |
| MW         | 1111.3                                                                     |
| Purity     | ≥97% (TLC), ≥92% (HPLC)                                                    |
| Solubility | Soluble in 100% ethanol (50mg/ml), acetic acid, chloroform, DMSO or ether. |
| Source     | Isolated from <i>Streptomyces fulvissimus</i>                              |

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-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)