

# Antibiotic A-23187 (Calcimycin)

Calcium ionophore

Antibiotic possessing weak *in vitro* antimicrobial activity against gram positive bacteria and fungi. A highly selective divalent cation ionophore, commonly used to increase intracellular  $\text{Ca}^{2+}$  levels in intact cells.

Citations: 12

[View Online »](#)

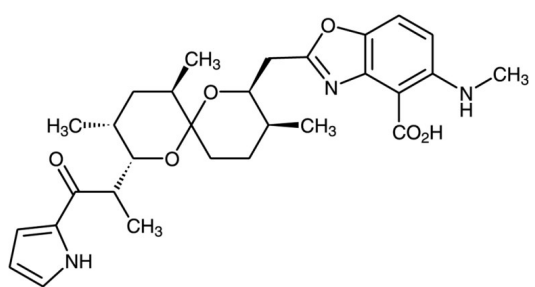
Ordering Information

[Order Online »](#)

BML-CA100-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. In aqueous systems product may aggregate over time. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
<b>Handling</b>	Protect from light.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Appearance</b>	Off-white solid.
<b>CAS</b>	52665-69-7
<b>Formula</b>	$C_{29}H_{37}N_3O_6$
<b>MW</b>	523.6
<b>Purity</b>	≥99% (TLC)
<b>Solubility</b>	Soluble in DMSO an, 100% ethanol or methylene chloride (10mg/ml); slightly soluble in water. Solubilize in DMSO or ethanol prior to preparing aqueous solutions of very low concentrations.
<b>Source</b>	Isolated from <i>Streptomyces chartreusis</i>



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-  
be@enzolifesciences.com](mailto:info-be@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)