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 Ca²⁺ levels in intact cells.

Citations: 12

View Online »

Ordering Information

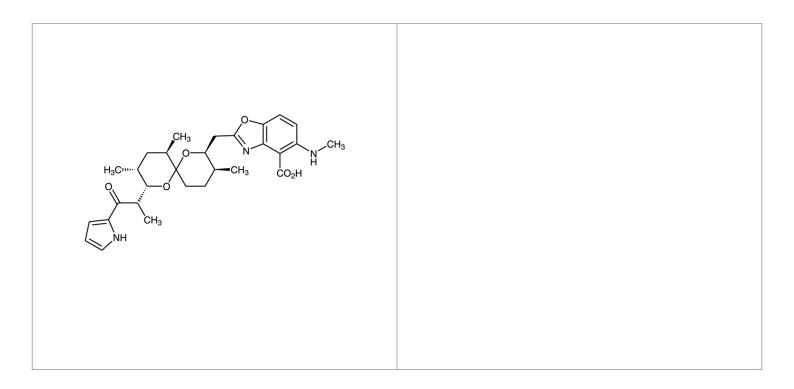
Order Online »

BML-CA100-0010

10mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability 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.

Handling Protect from light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Off-white solid.

CAS 52665-69-7

Formula $C_{29}H_{37}N_3O_6$

MW 523.6

Purity ≥99% (TLC)

Solubility 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.

Source Isolated from Streptomyces chartreusis

Last modified: May 29, 2024



European Sales Office