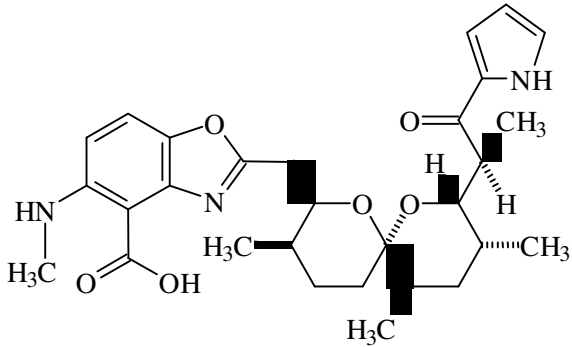


(also known as Calcimycin or Calcium Ionophore III)

Instruction Manual

Catalog Number	PK-CA707-59001
Description	A-23187 is a calcium ionophore that rapidly equilibrates intracellular and extracellular calcium concentrations. It is commonly used for in situ calibrations of fluorescent calcium indicators. The ionophore allows Mn^{2+} to enter the cells and quench intracellular fluorescence of calcium indicators.
Quantity	1 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	523.6 Da; $C_{29}H_{37}N_3O_6$
Purity	$\geq 98\%$ by TLC
Appearance / Formulation / Solubility	White solid; soluble in DMSO or EtOH.
Storage & Stability	Store at $-20^{\circ}C$ and protect from light, especially when in solution.
Applications	Calcium ionophor
References	<ol style="list-style-type: none"> 1. Suzuki, K. et al. Anal. Chem. 61, 382(1989) 2. Caron-Leslie, L.A., et al. FASEB J. 8, 639(1994) 3. Merck Index 12, 1678 4. Saunders, M.J., et al. Science 217, 943(1982)
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.