

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

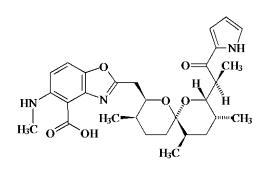
www.biotium.com

## **PRODUCT AND SAFETY DATA SHEET**

**PRODUCT NAME:** 

A-23187 Free Acid (also known as Calcimycin or Calcium Ionophore III)

**CATALOGUE #:** 59001



White solid $\geq$ 98% by TLC Soluble in DMSO or EtOH
Store at 4 °C and protect from light, especially when in solution.
A-23187 is a calcium ionophore that rapidly equilibrates intracellular and extracellular calcium concentrations. It is commonly used for <i>in situ</i> calibrations of fluorescent calcium indicators. The ionophore allows Mn <sup>2+</sup> to enter the cells and quench intracellular fluorescence of calcium indicators. Ref: 1) Suzuki, K. et al. <i>Anal. Chem.</i> <b>61</b> , 382 (1989); 2) Caron-Leslie, L.A., et al. <i>FASEB J.</i> <b>8</b> , 639 (1994); 3) <i>Merck Index</i> <b>12</b> , 1678; 4) Saunders, M.J., et al. <i>Science</i> <b>217</b> , 943 (1982).
DM4676000
LD50: 9.2 mg/kg, intraper, mouse; LD50: 4.5 mg/kg, intrav, mouse.
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.

**Disclaimer:** *Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.*