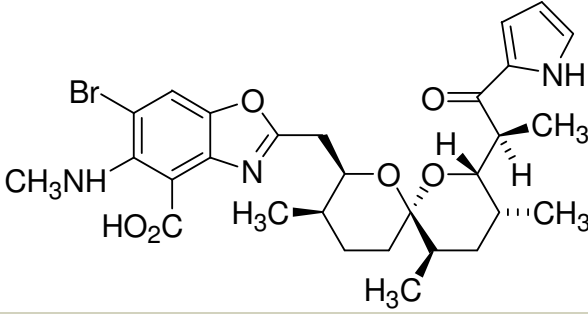


4-Bromo A-23187 free acid

For *in situ* calibration of fluorescent calcium indicators

Instruction Manual

| | |
|---------------------------------------|---|
| Catalog Number | PK-CA707-59006 |
| Description | Similar to A-23187, 4-Bromo A-23187 is used for <i>in situ</i> calibrations of fluorescent Ca^{2+} indicators to equilibrate intracellular and extracellular calcium concentrations and to allow Mn^{2+} to enter cells to quench indicator fluorescence. However, since 4-bromo A-23187 is nonfluorescent, it is preferred over the blue fluorescent A-23187 for calibrations of UV-excited Ca^{2+} indicators such as Indo-1 and Fura-2 in order to keep background fluorescence to a minimum. |
| Quantity | 1 mg |
| Excitation / Emission Maxima | NA |
| Molecular Structure |  |
| Molecular Weight / Molecular Formula | 602.52 Da; $\text{C}_{29}\text{H}_{36}\text{BrN}_3\text{O}_6$ |
| Purity | >98% (as determined by TLC) |
| Appearance / Formulation / Solubility | White solid; soluble in DMSO or EtOH. |
| Storage & Stability | Store at 4°C. Protect from light. |
| Applications | <i>In situ</i> calibrator for fluorescent Ca^{2+} indicators |
| References | Anal. Chem. 146, 349 (1985) |
| 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.