A-23187 Free Acid

A-23187, also known as Calcimycin or Calcium Ionophore III, is a calcium ionophore that rapidly equilibrates intracellular and extracellular calcium concentrations. It is commonly used for *in situ* calibrations of fluorescent calcium indicators.



Product attributes

CAS number

52665-69-7

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

A-23187, also known as Calcimycin or Calcium Ionophore III, 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 Mn² to enter the cells and quench intracellular fluorescence of calcium indicators. Also see 4-bromo A-23187 (59006). Because 4-bromo A-23187 is nonfluorescent, it is preferred over the blue fluorescent A-23187 for calibrations of UV-excited Ca² indicators such as Indo-1 and Fura-2 in order to keep background fluorescence to a minimum.

- White solid soluble in DMSO
- Store at 4°C and protect from light
- C29H37N3O6
- MW: 523.6
- [52665-69-7]

$$\begin{array}{c|c} CH_3NH & O & O & NH \\ HO_2C & H_3C & H_3C & CH_3 \end{array}$$

References

- 1. Anal Chem 61, 382 (1989).
- 2. FASEB J 8, 639 (1994).
- 3. Merck Index 12, 1678.
- 4. Science 217 (4563), 943 (1982).

This datasheet was generated on November 9, 2025 at 08:41:32 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/a-23187-free-acid-calcimycin-or-calcium-ionophore-iii/