

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.



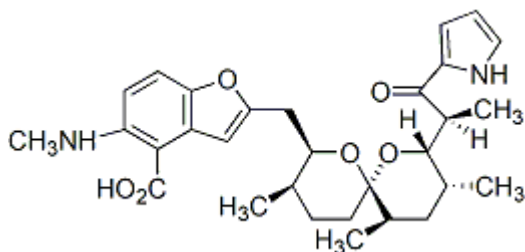
Product attributes

CAS number	52665-69-7
------------	------------

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 Mn^{2+} 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 Ca^{2+} 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
- $C_{29}H_{37}N_3O_6$
- MW: 523.6
- [52665-69-7]



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 June 8, 2026 at 04:59:03 AM. 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/>