

4-Bromo A-23187, Free Acid

Similar to A-23187, 4-Bromo A-23187 is used for in situ 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.



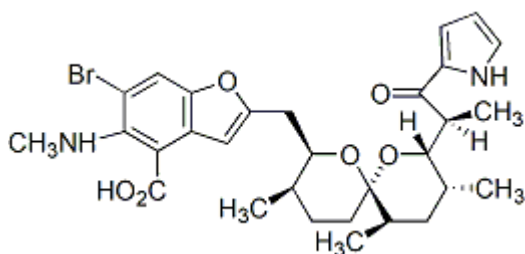
Product attributes

CAS number	76455-48-6
------------	------------

Product Description

Similar to A-23187, 4-Bromo A-23187 is used for in situ 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.

- White solid soluble in DMSO or EtOH
- Store at 4 °C
- $\text{C}_{29}\text{H}_{36}\text{BrN}_3\text{O}_6$
- MW: 602.52
- [76455-48-6]



References

- 1) Anal Chem 146, 349 (1985).

This datasheet was generated on June 25, 2026 at 09:25:53 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/4-bromo-a-23187-free-acid/>