Ionomycin, Calcium Salt

lonomycin is commonly used to modify intracellular calcium concentrations and to calibrate fluorescent calcium indicators. It is also known for its antiproliferative effects.



Product attributes

CAS number

56092-82-1

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

Product Description

lonomycin is commonly used to modify intracellular calcium concentrations and to calibrate fluorescent calcium indicators. It is also known for its antiproliferative effects.

- Yellowish oily solid soluble in DMSO or EtOH
- Store desiccated at ≤ 4°C
- C₄₁H₇₀O₉Ca
- MW: 747.1
- [56092-82-1]

References

- 1) J Biol Chem 275, 7071 (2000).
- 2) J Immunol 155, 3297 (1995).
- 3) J Urol 162, 916 (1999).

This datasheet was generated on November 3, 2025 at 03:53:18 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/ionomycincalcium-salt/