

## Ionomycin, Calcium Salt

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



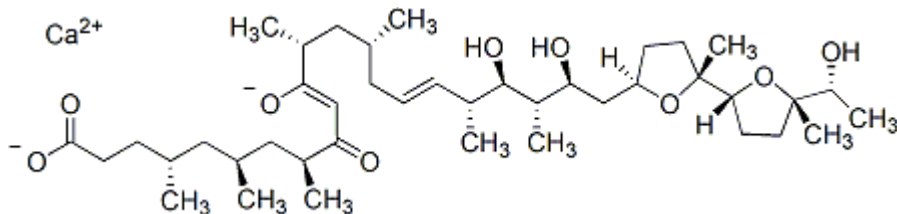
### Product attributes

CAS number	56092-82-1
------------	------------

## Product Description

Ionomycin 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  $\leq 4^{\circ}\text{C}$
- $\text{C}_{41}\text{H}_{70}\text{O}_9\text{Ca}$
- 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 June 7, 2026 at 10:11:47 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/ionomycincalcium-salt/>