

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

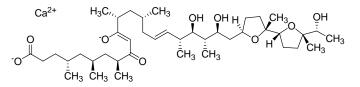
PRODUCT NAME:

Ionomycin, calcium salt

CATALOG #

MOLECULAR INFORMATION: C₄₁H₇₀O₉Ca Mwt: 747.1 CAS: [56092-82-1]

59007



PROPERTIES:

Color & Form	Off-white to yellowish solid
Solubility	Soluble in DMSO or EtOH

STORAGE AND Store desiccated at \leq 4 °C. **HANDLING:**

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

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

TOXICITY: LD₅₀: 650 mg/kg, oral, mouse; 12 mg/kg, intraperitoneal, mouse.

FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: *Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.*