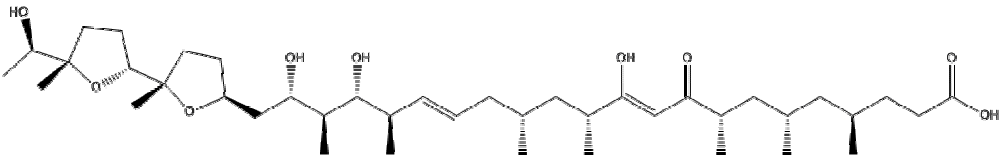


## Instruction Manual

Catalog Number	PK-CA577-1565-5
Description	Antibiotic that acts as a potent and selective calcium ionophore; more effective than A23187. Ionomycin binds $\text{Ca}^{2+}$ in the 7.0-9.5 pH range with the resulting complex exhibiting intense UV absorption. Ionomycin induces apoptosis in immature B cell lines and in cultured embryonic rat cortical neurons. More mature B cell lines do not undergo apoptosis, but are arrested in the G1 phase of the cell cycle.
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	709.02 Da; $\text{C}_{41}\text{H}_{72}\text{O}_9$
Chemical Structure	
Purity	>97%
Appearance / Formulation / Solubility	Waxy solid or off-white to light yellow powder. Soluble in DMSO.
Storage & Stability	-20°C in the dark. Protect from light and moisture. Do not expose to sunlight or fluorescent lights for extended periods.
Applications	see Description
References	NA
Caution	Tumor promoter. Do not ingest, swallow, or inhale. Do not get in eyes or on skin. Care should be taken to prevent contact through all routes of exposure.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.