

Instruction Manual

Catalog Number	PK-CA577-1558-1
Description	Potent, cell permeable IP3-independent intracellular calcium releaser; Inhibits microsomal Ca ²⁺ -ATPase; Stimulates arachidonic acid metabolism in macrophages.
Quantity	1 mg
Sequence / Molecular Weight / Molecular Formula	650,76 Da; C ₃₄ H ₅₀ O ₁₂
Purity	>99%
Appearance / Formulation / Solubility	Powder. Soluble in DMSO or 100% ethanol.
Storage & Stability	Store at -20°C.
Applications	see Description
References	NA
Caution	Irritant to skin and mucous membranes. Possible tumor promoter. Wear gloves and mask when using this compound. Care should be taken to prevent contact through all routes of exposure.

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