

Instruction Manual

Catalog Number	PK-CA577-1544-5
Description	This is the most commonly used phorbol ester. It binds to and activates protein kinase C, causes an extremely wide range of effects in cells and tissues and is a very potent mouse skin tumor promoter. Inhibits apoptosis induced by the Fas antigen, but induced apoptosis in HL-60 promyelocytic leukemia cells.
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	616,8 Da; C ₃₅ H ₅₅ O ₅
Purity	>99%
Appearance / Formulation / Solubility	Soluble in DMSO or ethanol.
Storage & Stability	Store at -20°C. Protect from light!
Applications	see Description
References	NA
Caution	Highly irritant 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. Toxicity information is not available at this time.

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