

Instruction Manual

Catalog Number	PK-CA577-1560-5
Description	Brefeldin A has been shown to induce apoptosis in human tumor cells through caspase activation. It has also been shown to block translocation of proteins from the endoplasmic reticulum to the Golgi apparatus.
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	280,4 Da; C ₁₅ H ₂₄ O ₄
Purity	>99%
Appearance / Formulation / Solubility	powder. Soluble in DMSO and ethanol.
Storage & Stability	Store at -20°C in the dark.
Applications	see Description
References	NA
Caution	Toxic. 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. Detailed toxicity information is not available at this time.

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