

## Instruction Manual

Catalog Number	PK-CA577-1519
Description	Inhibitor of microtubules by specific binding to tubulin. Antimitotic tool used in cell division studies; Antibiotic; anti-inflammatory; therapeutic agent. Induce apoptosis in PC-12 cells and cerebellar granule cells.
Quantity	1 g
Sequence / Molecular Weight / Molecular Formula	399.4 Da; C <sub>22</sub> H <sub>25</sub> NO <sub>5</sub>
Purity	>99%
Appearance / Formulation / Solubility	White solid. Soluble in DMSO and water
Storage & Stability	Store desiccated at -20°C in the dark.
Applications	see Description
References	1. Sherline, P., et al. (1974) Anal. Biochemistry 62:400. 2. Arai, T. and Okuyama, T. (1975) Anal. Biochemistry 69:443.
Caution	Potent neurotoxin!

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