

## Instruction Manual

Catalog Number	PK-CA577-1524-10
Description	Potent anti-cancer agent whose potential target site may be mitochondrial cytochrome c oxidase. Inhibits RNA and DNA synthesis. Also inhibits eukaryotic topoisomerase I & II. Induces DNA single strand breaks and apoptosis in HeLaS3 tumor cells.
Quantity	10 mg
Sequence / Molecular Weight / Molecular Formula	564 Da; C <sub>25</sub> H <sub>29</sub> NO <sub>10</sub> .HCl
Purity	>98%
Appearance / Formulation / Solubility	Orange-Red Power. Soluble in water or methanol.
Storage & Stability	Store at -20°C in a dry environment.
Applications	see Description
References	1. Bose, R., et al. (1995) Cell 82:405. 2. Boland, M.P., et al. (1997) J. Biol. Chem. 272:12952.
Caution	NA

---

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

---