

## Instruction Manual

Catalog Number	PK-CA577-1042-1
Description	An anti-inflammatory glucocorticoid. Inhibits the expression of the inducible but not the constitutive nitric oxide synthase in vascular endothelial cells (IC50=5 nM). Enhances active cation transport in aortic smooth muscle cells by stimulating the Na <sup>+</sup> -K <sup>+</sup> pump. Has antiinflammatory and anti-rheumatic properties. Induces apoptosis in human thymocytes. In general, 500-1000 nM of dexamethasone is sufficient to induce apoptosis following a 6-hour incubation at 37°C (1, 2).
Quantity	1 ml
Sequence / Molecular Weight / Molecular Formula	392,5 Da; C <sub>22</sub> H <sub>29</sub> FO <sub>5</sub>
Purity	>99% by TLC
Appearance / Formulation / Solubility	10 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 3 months under proper storage conditions.
Applications	We recommend using 1000X dilutions for inducing apoptosis in Jurkat cells. However, the optimal doses may vary for different cells and culture conditions.
References	1. Martin, S.J., et al. (1995) J. Exp. Med. 182:1545-1556. 2. Fearnhead, H.O., et al. (1994) Biochem. Pharmacol. 48:1073-1079.
Caution	NA

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