

## Instruction Manual

Catalog Number	PK-CA577-1568-50
Description	Immunosuppressant. Selectively blocks signaling that leads to p70 S6 kinase activation (IC <sub>50</sub> = 50 pM). Inhibits lymphokine-induced cell proliferation at the G1 phase. Induce apoptosis in a murine B cell line. Rapamycin has been used in vivo studies by i.p., i.m., and i.v. administrations.
Quantity	50 mg
Sequence / Molecular Weight / Molecular Formula	914,19 Da; C <sub>51</sub> H <sub>59</sub> NO <sub>13</sub>
Purity	>99%
Appearance / Formulation / Solubility	White to off-white solid. Soluble in DMSO and ethanol.
Storage & Stability	-20°C in the dark. Do not expose to sunlight or fluorescent lights for extended periods.
Applications	see Description
References	NA
Caution	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.

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