

Rapamycin (Sirolimus)

Catalog #: 27062

Size: 50 mg

CAS Registry #: 53123-88-9

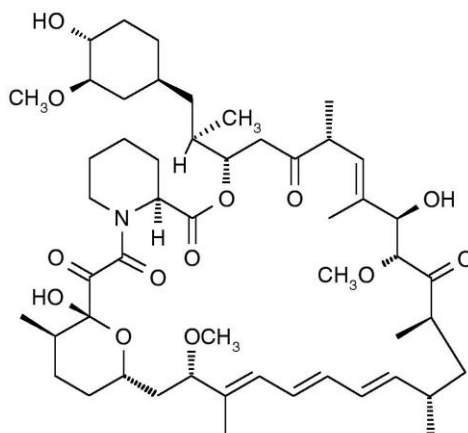
Purity: ≥99%

Chemical Formula: C₅₁H₇₉NO₁₃

Molecular Weight: 914.17 Da

Lot #: 230306

Structure:



Description: Immunosuppressant, related to FK-506, but without calcineurin inhibitory activity even when complexed to FK-506 binding protein. Selectively blocks signaling that leads to p70 S6 kinase activation (IC₅₀ = 50 pM).

Appearance: White to off-white crystalline solid, granular or powder

Solubility: Soluble in DMSO at 200 mg/ml; soluble in ethanol at 50 mg/ml; very poorly soluble in water; maximum solubility in plain water is estimated to be about 5-20 μM; buffers, serum, or other additives may increase or decrease the aqueous solubility.

Storage/Stability: Store at or below -20°C. Solid form is stable at least 12 months from date of receipt, when stored as directed. Do not store aqueous solutions for more than one day. *Avoid freeze/thaw cycles.*

References:

1. Terada, N., *et al.*, *J. Biol. Chem.* 1993; **268**: 12062-12068.