FK506

Immunosuppressant

Immunosuppressant that blocks T cell proliferation *in vitro* by inhibiting the generation of several lymphokines, especially IL-2. Shown to inhibit the activity of FK506 binding protein and to inhibit the Ca^{2+} -dependent phosphatase and protein phosphatase 2B (PP2B; calcineurin). Activates NF- κ B through phosphorylation and degradation of $I\kappa$ B α .

Citations: 27

View Online »

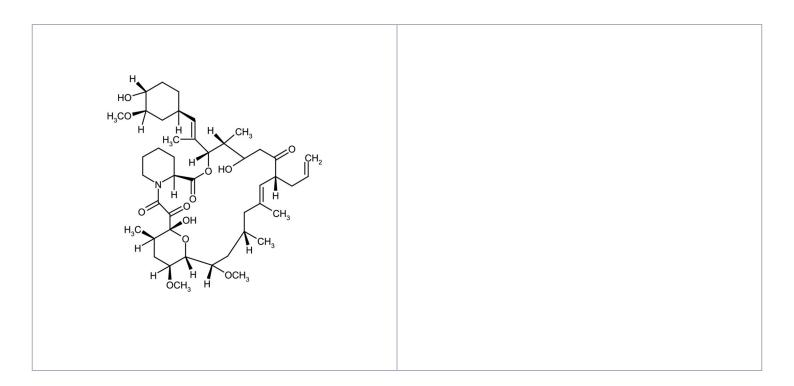
Ordering Information

Order Online »

ALX-380-008-M005	5mg
ALX-380-008-M025	25mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tacrolimus

Appearance White to off-white crystalline powder.

CAS 104987-11-3

Couple Target FKBP, Serine/threonine-protein phosphatase

Couple Type Inhibitor, Ligand

Formula $C_{44}H_{69}NO_{12}$

Identity FT-IR

14: 9025 ΜI

MW 804.0

Purity ≥98% (HPLC)

RTECS KD4200000

Solubility Soluble in DMSO, 100% ethanol, methanol, ethyl acetate

or acetone; insoluble in water.



Phone: 800.942.0430 infousa@enzolifesciences.com

ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 eu@enzolifesciences.com

European Sales Office

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Isolated from Streptomyces tsukubaensi. Phone: +33 472 440 655 fr@enzolifesciences.com

Phone: +49 7621 5500 526 de@enzolifesciences.com

0845 601 1488 Phone: +44 1392 825900 uk@enzolifesciences.com

Last modified: May 29, 2024