Anisomycin

Activator of MAPK, Jnk and p38

Antibiotic. Activates JNK/SAPKs, p38 and MAP kinases but not ERKs leading to the phosphorylation of S6, histone H3 and HMG-14 and the induction of the immediate-early genes c-fos and c-jun. Induces apoptosis in the human monoblastoid cell line. Used at 300 ng/ml. Inhibits peptidyl transferase at higher concentrations. Used in the eradication of bean mildew. Inhibits other pathogenic fungi in plants.

Citations: 18

View Online »

Ordering Information

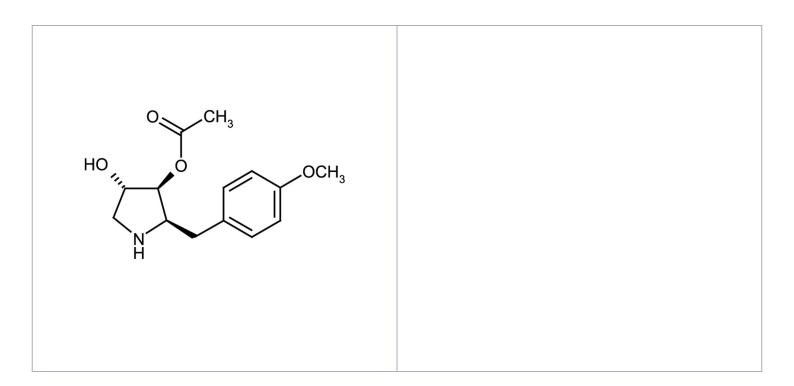
Order Online »

BML-ST102-0050

50mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up

to 3 months at -20°C.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-(p-Methoxybenzyl)-3,4-pyrrolidinediol-3-acetate

Appearance White solid.

CAS 22862-76-6

Couple Target Jnk, MAPK, p38

Couple Type Activator

Formula $C_{14}H_{19}NO_4$

MI 14: 670

MW 265.3

Purity ≥97% (TLC)

RTECS BZ9800000

Soluble in DMSO (25mg/ml).

Source Isolated from Streptomyces griseolus.