S14-95

STAT1α inhibitor

Potent inhibitor of cytokine-induced activation of STAT1 α leading to the inhibition of inducible expression of proinflammatory enzymes (COX-2, iNOS (NOS II)) and cytokines (TNF- α).

Citations: 4

View Online »

Ordering Information

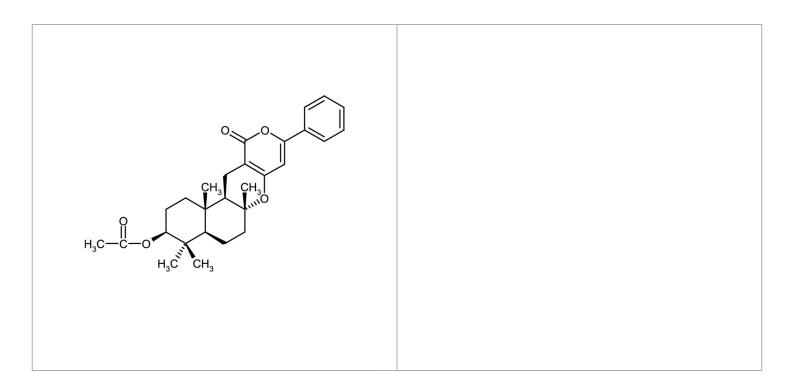
Order Online »

ALX-350-299-MC05

0.5mg

Manuals, SDS & CofA

View Online »



Handling & Storage

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

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Yellow powder.

Couple Target STAT

Couple Type Inhibitor

Formula $C_{28}H_{34}O_5$

Identity Determined by MS.

MW 450.6

Purity ≥95% (HPLC)

Solubility Soluble in methanol.

Source Isolated from *Penicillium* sp.

Last modified: May 29, 2024