

S14-95

STAT1 α 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 »](#)

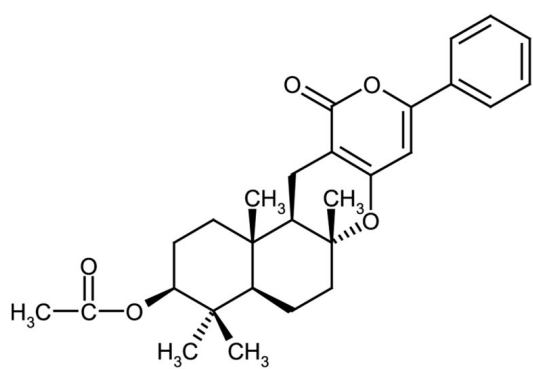
Ordering Information

[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 <i>Penicillium</i> sp.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)