

Galiellalactone

STAT3 inhibitor

Originally isolated from *Galiella rufa* as a plant growth regulator. Inhibits IL-6 mediated signal transduction by blocking the binding of the activated STAT3 dimers to their DNA binding sites without inhibiting the tyrosine and serine phosphorylation site of the STAT3 transcription factor.

Citations: 7

[View Online »](#)

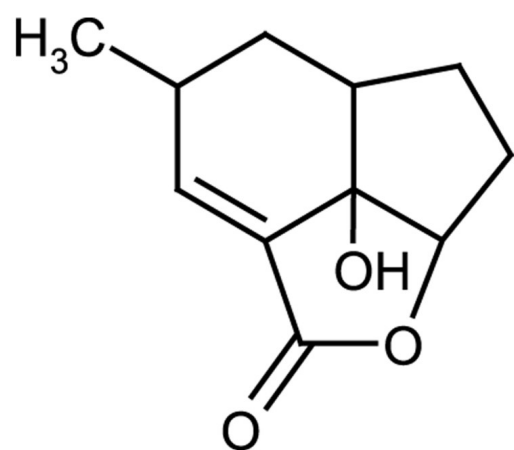
Ordering Information

[Order Online »](#)

ALX-350-336-MC05	0.5mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to light tan solid.

CAS 133613-71-5

Couple Target STAT

Couple Type Inhibitor

Formula $C_{11}H_{14}O_3$

Formulation Lyophilized.

MW 194.2

Purity ≥95% (HPLC)

Solubility Soluble in 100% ethanol, methanol, dimethyl formamide or DMSO.

Source Isolated from an unidentified fungus, MST-FP1889.

Last modified: May 29, 2024



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)