

Rocaglamide

NF-κB inhibitor

Immunosuppressant. Potent inhibitor of NF-κB activation in T cells, with an almost complete inhibition at 200nM. Suppresses cytokine production (IFN-γ, TNF-α, IL-2 and IL-4) and inhibits NF-AT in peripheral blood T cells at concentrations that do not impair NF-κB and AP-1 activities. In contrast to the immunosuppressant cyclosporin A (Prod. No. BML-A195), rocaglamide does not inhibit calcineurin phosphatase activity. Induces apoptosis.

Citations: 17

[View Online »](#)

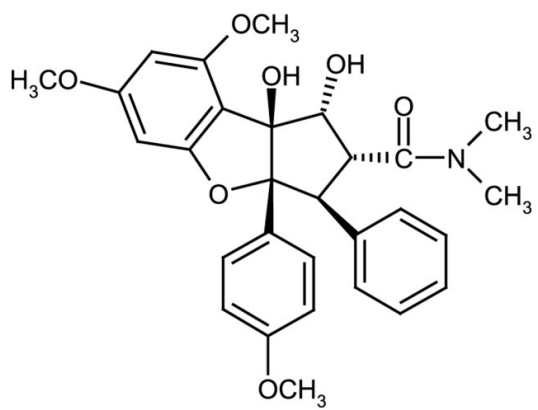
Ordering Information

[Order Online »](#)

ALX-350-121-C100	100μg
------------------	-------

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	Rocaglamide A
Appearance	White to off-white sticky solid.
CAS	84573-16-0
Couple Target	NF-kappaB
Couple Type	Inhibitor
Formula	$C_{29}H_{31}NO_7$
MW	505.6
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO, 100% ethanol or methanol.
Source	Synthetic

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-be@enzolifesciences.com)

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