

Ratjadone A

(synthetic)

Represents a class of natural compounds, which inhibit proliferation in eukaryotes by blocking nuclear export. As potent as leptomycin B (Prod. No. ALX-380-100) and specific for G1/S checkpoint. Cytotoxic secondary metabolite (IC_{50} =50pg/ml in mouse cell line L929) that arrests tumor cells in the G1 phase at remarkably low concentrations (50pg/ml in HeLa cell line KB3.1). Inhibits the binding between the nuclear export signal (NES) of proteins and the chromosome maintenance region protein (CRM1). Anticancer compound. Belongs to the family of orphan ligands which include polyketides like leptomycin B, callistatin A and other related compounds.

Citations: 6

[View Online »](#)

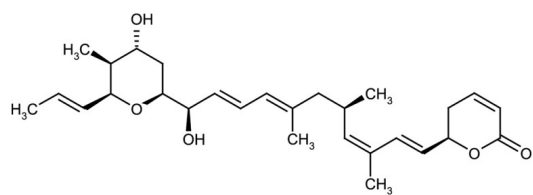
Ordering Information

[Order Online »](#)

ALX-270-346-C002	2µg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Sensitive to acids and bases. Unstable in acidic buffer systems (pH<6). Stable in substance as oil and in solution in methanol at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Clear liquid.
CAS	163564-92-9
Formula	$C_{28}H_{40}O_5$
Formulation	Solution in methanol.
Identity	Determined by MS, 1H- and 13C-NMR.
MW	456.6
Purity	≥95% (HPLC)
Solubility	Soluble in methanol (20mg/ml) or aqueous buffers (<100µM).
Source	Synthetic.



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)