

Avarol

Cytostatic agent

Cytostatic agent which has potent antileukemic activity both *in vitro* and *in vivo* (mice). Also displays antibacterial and antifungal activities against a limited range of microorganisms. Inhibits HIV-1 reverse transcriptase.

Citations: 20

[View Online »](#)

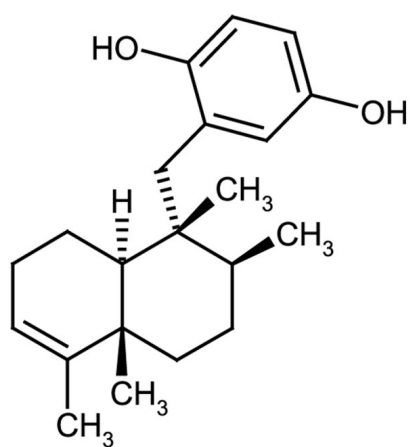
Ordering Information

[Order Online »](#)

ALX-350-319-M001	1mg
------------------	-----

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	NSC306951
Appearance	Brown solid.
CAS	55303-98-5
Couple Target	Reverse transcriptase
Couple Type	Inhibitor
Formula	$C_{21}H_{30}O_2$
Identity	Determined by 1H -NMR, ^{13}C -NMR and MS.
MW	314.5
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO or 100% ethanol.
Source	Isolated from marine sponge <i>Dysidea avara</i> .

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)