

# Avarone

## 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.

Citations: 13

[View Online »](#)

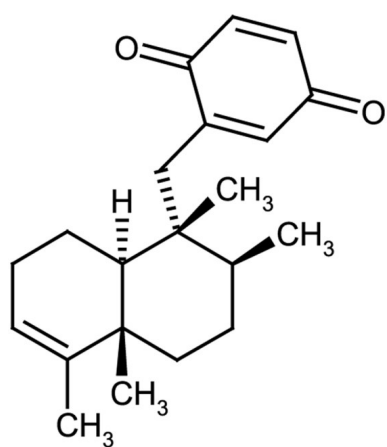
## Ordering Information

[Order Online »](#)

ALX-350-321-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

**Appearance** Brown oil.

**CAS** 55303-99-6

**Formula**  $C_{21}H_{28}O_2$

**Identity** Determined by <sup>1</sup>H-NMR, <sup>13</sup>C-NMR and MS.

**MW** 312.5

**Purity** ≥97% (HPLC)

**Solubility** Soluble in DMSO or 100% ethanol.

**Source** Isolated from marine sponge *Dysidea avara*.

**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)