

# Gilvocarcin M

Antitumor antibiotic

Antibiotic. Has antibacterial, antifungal, antiviral and antitumor activities. Interacts with DNA. Less active than gilvocarcin V.

Citations: 3

[View Online »](#)

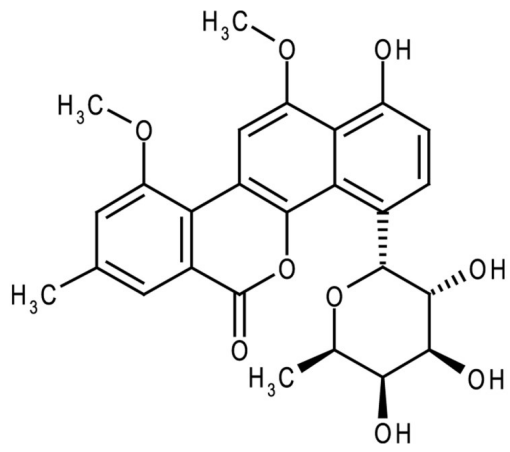
Ordering Information

[Order Online »](#)

ALX-380-228-MC05	0.5mg
------------------	-------

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	Anandimycin B, Toromycin B, Antibiotic 1072A
Appearance	Yellow solid.
CAS	77879-89-1
Formula	$C_{26}H_{26}O_9$
MW	482.5
Purity	≥95% (HPLC)
RTECS	DI2250000
Solubility	Soluble in dimethyl formamide or DMSO; slightly soluble in methanol or 100% ethanol.
Source	Isolated from <i>Streptomyces</i> sp. MST-AS5353.

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)