

Gilvocarcin M

Antitumor antibiotic

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

Citations: 3

[View Online »](#)

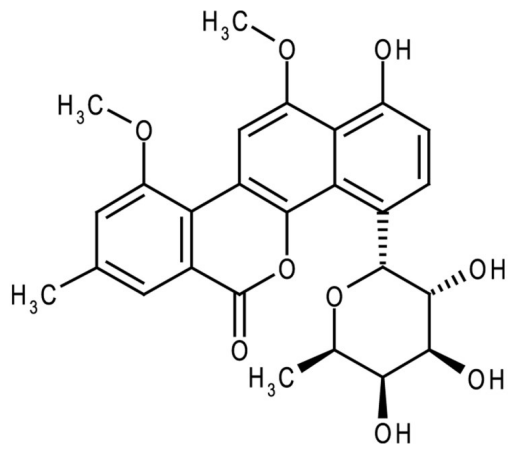
Ordering Information

[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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com