

# Meleagrín

Antibiotic

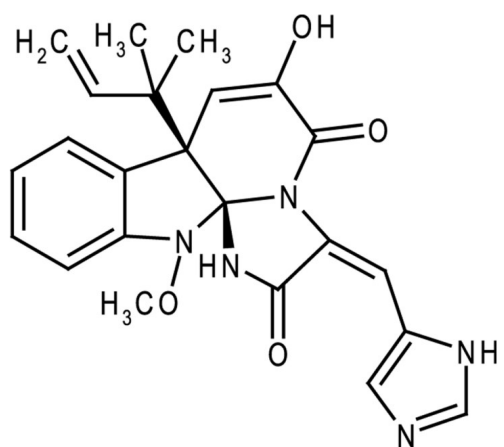
Antibiotic.

Citations: 3

[View Online »](#)

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 | Yellow oil.                                    |
| CAS        | 71751-77-4                                     |
| Formula    | $C_{23}H_{23}N_5O_4$                           |
| Identity   | Determined by 1H-NMR and MS.                   |
| MW         | 433.5  |
| Purity     | ≥98% (HPLC)                                    |
| Solubility | Soluble in 100% ethanol or DMSO.               |
| Source     | Isolated from <i>Penicillium meleagrinum</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-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)