

Lycorine . HCl

Protein synthesis inhibitor

Lycorine is a eukaryotic termination inhibitor.

Citations: 2

[View Online »](#)

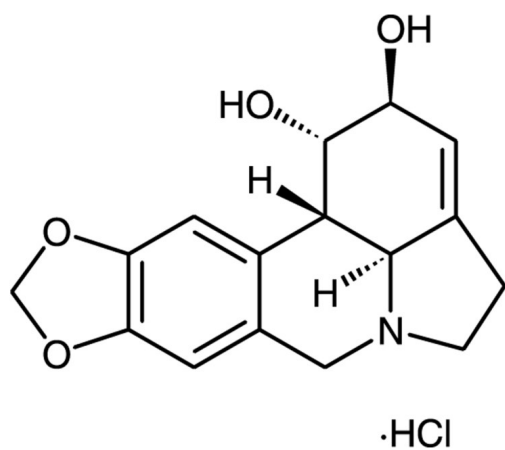
Ordering Information

[Order Online »](#)

BML-GR313-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white powder.
CAS	2188-68-3
Formula	$C_{16}H_{17}NO_4 \cdot HCl$
Identity	Determined by NMR.
MW	323.8
Purity	≥97% (Assay)
Solubility	Soluble in water (>15mg/ml).

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