

Gliotoxin

Proteasome inhibitor

Immunomodulating mycotoxin which acts by blocking membrane thiol groups. Inhibits the chymotrypsin-like activity of the 20S proteasome in a non-competitive manner. Causes apoptotic cell death in macrophages and thymocytes. Induces Ca^{2+} release from intact rat liver mitochondria. Inhibits the activation of transcription factor NF- κ B in response to a variety of stimuli in T and B cells. Displays anti-inflammatory activity *in vivo*. Antioxidant. Inhibitor of farnesyltransferase and geranylgeranyltransferase I. A negative control analog is available (Prod. No. BML-L122). Displays anti-tumor activity against breast cancer *in vivo*.

Citations: 13

[View Online »](#)

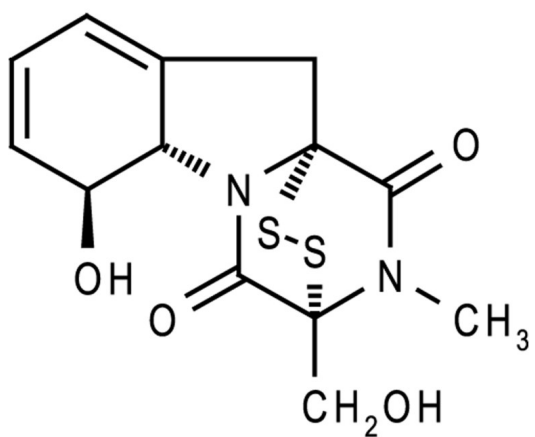
Ordering Information

[Order Online »](#)

BML-PI129-0002	2mg
BML-PI129-0010	10mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	[3R-(3 α , 5 $\alpha\beta$, 6 β , 10 $\alpha\alpha$)]-2,3,5a,6-Tetrahydro-6-hydroxy-3-(hydroxymethyl)-2-methyl-10H-3,10a-epi-dithiopyrazino[1,2-a]indole-1,4-dione
Appearance	White to yellow crystalline solid.
CAS	67-99-2
Couple Target	Proteasome 20S
Couple Type	Inhibitor
Formula	C ₁₃ H ₁₄ N ₂ O ₄ S ₂
MI	14: 4441
MW	326.4
Purity	≥97% (HPLC, TLC)
RTECS	KB4725000
Solubility	Soluble in DMSO or methanol (10mg/ml).
Source	Isolated from <i>Gliocladium fimbriatum</i>



ENZO LIFE SCIENCES,
Last modified: **May 29, 2024**
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