

Mitomycin C

DNA synthesis inhibitor

Antitumor antibiotic. DNA crosslinking and damaging agent. Inhibitor of DNA synthesis, nuclear division and cancer cells. DNA intercalating agent. Induces apoptosis in a variety of cells via a p53-dependent pathway.

Citations: 28

[View Online »](#)

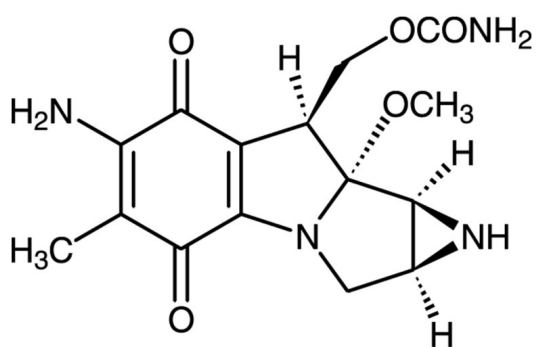
Ordering Information

[Order Online »](#)

| | |
|----------------|------|
| BML-GR311-0002 | 2mg |
| BML-GR311-0010 | 10mg |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|-------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -20°C. Stock solutions in water or ethanol are stable for up to 1 week at -20°C. Stock solutions in DMSO are stable for up to 1 month at -20°C. |
| Handling | Protect from light. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|---------------|---|
| Appearance | Blue-violet solid. |
| CAS | 50-07-7 |
| Couple Target | DNA |
| Couple Type | Ligand |
| Formula | $C_{15}H_{18}N_4O_5$ |
| MI | 14: 6215 |
| MW | 334.3 |
| Purity | ≥95% (HPLC) |
| RTECS | CN0700000 |
| Solubility | Soluble in water (1mg/ml, warm), 100% ethanol (1mg/ml, warm), DMSO (15mg/ml) or methanol. |

Technical Info / Product Note: Product is not sterile.

Notes



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-fr@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