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 »

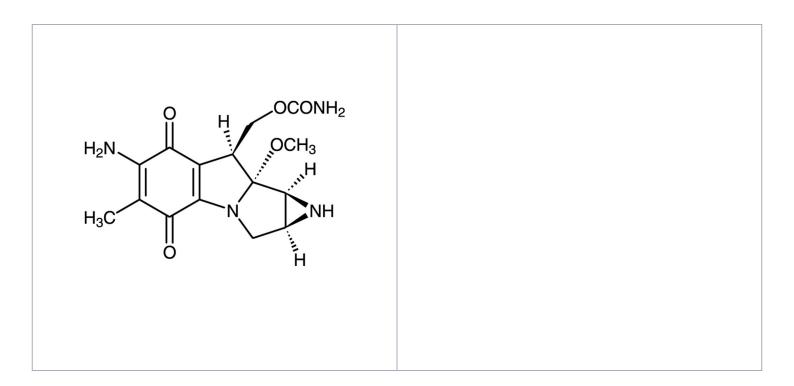
Ordering Information

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

Soluble in water (1mg/ml, warm), 100% ethanol (1mg/ml,

warm), DMSO (15mg/ml) or methanol.

Technical Info / Product NotesNote: Product is not sterile.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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