

Bleomycin sulfate

Antitumor antibiotic

Bleomycin consists in a complex of 11 glycopeptide anti-cancer antibiotics originally isolated from *Streptomyces verticillus*. The dominant components of the complex are bleomycin A2 and B2, which typically represent >90% of the total weight. Bleomycin has found clinical application in the treatment of a range of tumors. Bleomycin acts by intercalation of DNA and RNA. In the presence of oxygen and metal ions (e.g. copper and iron), Bleomycin forms a pseudo-enzyme that induces DNA cleavage.

Citations: 38

[View Online »](#)

Ordering Information

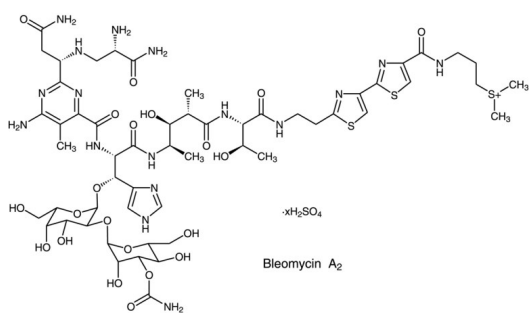
[Order Online »](#)

BML-AP302-0010	10mg
BML-AP302-0050	50mg

Manuals, SDS & CofA

[View Online »](#)

- Antibiotic with anti-cancer and apoptotic properties
- Radiomimetic DNA-cleaving agent that produces double and single DNA strand breaks through an oxygen-radical-dependent mechanism
- Highly cited



Bleomycin A₂



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, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Blenoxane
Appearance	White crystalline solid.
CAS	9041-93-4
Formula	$C_{55}H_{85}N_{17}O_{25}S_4$
MI	14: 1318
MW	1512.6
Purity	≥98% (HPLC)
RTECS	EC5991990
Solubility	Soluble in 100% methanol or water (20mg/ml).
Source	Isolated from <i>Streptomyces</i> sp.

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-
fr@enzolifesciences.com](mailto:info-fr@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)