

Bleomycin A5 . hydrochloride

Glycopeptide, induces DNA strand breaks.

Bleomycin A5 (pingyangmycin) is an antitumor antibiotic. Induces apoptosis.

Ordering Information

[Order Online »](#)

LKT-B4517-M010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Pingyangmycin hydrochloride

CAS 55658-47-4

Couple Type Inhibitor

Formula $C_{57}H_{89}N_{19}O_{21}S_2 \cdot HCl$

MW 1477.02

Purity $\geq 94\%$

Solubility Freely soluble in water, slightly soluble in ethanol



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