

Chromomycin A3

Transcription inhibitor

Chromomycin A₃ is a GC-specific DNA ligand which inhibits transcription. It may be used as a fluorescent DNA stain. Antibiotic. Cell permeable fluorescent dye that can be used with Bisbenzamide H33258 to distinguish chromosomes by their total DNA concentration and their DNA base composition. Does not intercalate. Ex(max): 445nm; Em(max): 575nm

Citations: 4

[View Online »](#)

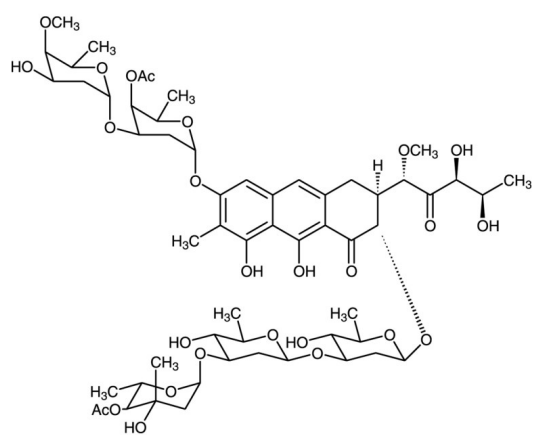
Ordering Information

[Order Online »](#)

BML-GR302-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow solid.
CAS	7059-24-7
Couple Target	DNA
Couple Type	Ligand
Formula	$C_{57}H_{82}O_{26}$
MW	1183.3
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol, DMSO or ethyl acetate (10mg/ml)
Source	Isolated from <i>Streptomyces griseus</i>



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