

G418 . sulfate

Antibiotic

G418 is a gentamycin-class aminoglycoside antibiotic with cytotoxic properties in both prokaryotic and eukaryotic cells. In a cell-free assay, G418 inhibits 80S ribosome and protein elongation during translation. In cells, G418 induces apoptosis and is widely employed in the selection of eukaryotic expression vectors containing the neomycin resistance gene neoR, in combination with either APH(2'') or APH(3').

This product is the disulfide salt form of G418 CAS# 49863-47-0.

Citations: 34

[View Online »](#)

Ordering Information

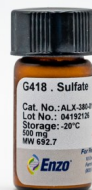
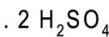
[Order Online »](#)

ALX-380-013-M500	500mg
ALX-380-013-G001	1g
ALX-380-013-G005	5g

Manuals, SDS & CofA

[View Online »](#)

- Used to select neoR-positive cell lines and plasmids
- Highly cited



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions are stable for 3 to 6 months when stored at -20°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Activity	Potency: ≥700 µg/mg
Alternative Name	Geneticin
Appearance	White to off-white powder.
CAS	108321-42-2
Formula	$C_{20}H_{40}N_4O_{10} \cdot 2H_2SO_4$
Identity	Identity determined by HPLC and IR.
MW	496.6 . 196.1
RTECS	CB9378500
Solubility	Soluble in water.
Source	Isolated from <i>Micromonospora rhodorangea</i> .

Technical Info / Product Notes	The effective killing concentration of this antibiotic will vary as to the cell type, media, growth conditions, and the cells metabolic rate and position in the cell cycle. Effective concentrations have been reported from 100µg/ml up to 5mg/ml or greater. When using G418 in a new cell system, a full dose curve is suggested and with each new lot of G418, several points on that curve should be retested, as the potency determined in a standard B. subtilis assay may not exactly correlate to your system.
--------------------------------	--



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)