

CELL CULTURE

SELECTIVE ANTIBIOTICS



Selection starts here

- ❖ Cell culture ready
- ❖ Endotoxin tested & rigorous quality control
- ❖ The highest quality at the best price

Discover the most extensive line of selective antibiotics you'll find anywhere. InvivoGen provides high-quality and affordable antibiotics for selection with unmatched purity levels. Manufactured in its state-of-the-art facility from its own proprietary strains, all InvivoGen selective antibiotics are stable, cell culture-tested and ready to use.

The selective antibiotics of choice

Blasticidin

G418 Sulfate

Hygromycin B Gold

Pleomycin

Puromycin

Zeocin™

WWW.INVIVOGEN.COM/SELECTIVE-ANTIBIOTICS

Selective Antibiotics

• Blasticidin

This nucleoside antibiotic specifically inhibits protein synthesis by interfering with the peptide-bond formation in the ribosomal machinery in both prokaryotes and eukaryotes. Resistance is conferred by the blasticidin resistance gene, bsr, derived from *Bacillus cereus*.

Purity: ≥95%

• G418 Sulfate

This aminoglycoside antibiotic blocks polypeptide synthesis by inhibiting the elongation step in both prokaryotes and eukaryotes. Resistance is conferred by the neo gene from Tn5.

Purity: ≥90%

• Hygromycin B Gold

This amino-glycoside antibiotic inhibits protein synthesis by interfering with translocation and causing mistranslations at the 70S ribosome. It is effective on most bacteria, fungi and higher eukaryotes. Resistance is conferred by the hph gene from *E. coli*.

Purity: ≥90%

• Phleomycin

A glycopeptide antibiotic of the bleomycin family, Phleomycin binds and intercalates DNA in most aerobic cells, destroying the integrity of the double helix. Phleomycin is active against most bacteria, filamentous fungi, yeast, plant and animal cells. Resistance is conferred by the Sh ble gene from *Streptoalloteichus hindustanus*.

• Puromycin

This aminonucleoside antibiotic inhibits peptidyl transfer on both prokaryotic and eukaryotic ribosomes. It inhibits the growth of Gram positive bacteria and various animal and insect cells. Resistance is conferred by the Pac gene.

Purity: ≥98%

• Zeocin™

A copper-chelated glyco-peptide antibiotic that causes cell death by intercalating into DNA and cleaving it. Effective on most aerobic cells, Zeocin™ is useful for selection in bacteria, eukaryotic microorganisms, plant and animal cells. Resistance is conferred by the Sh ble gene.

PRODUCT	WORKING CONC.	STABILITY	CONCENTRATION*	QUANTITY	CAT. CODE
Blasticidin	Cells: 1-10 µg/ml E. coli: 25-100 µg/ml	3 years at -20°C 2 years at 4°C 3 months at 20-25°C	10 mg/ml	50 mg (5 x 1 ml) 100 mg (10 x 1 ml) 500 mg (50 x 1 ml) 500 mg (50 ml bottle) 1 g (powder)	ant-bl-05 ant-bl-1 ant-bl-5 ant-bl-5b ant-bl-10p
G418 Sulfate	Cells: 400-1000 µg/ml	2 years at -20°C 1 year at 4°C 3 months at 20-25°C	100 mg/ml	1 g 5 g	ant-gn-1 ant-gn-5
Hygromycin B Gold	Cells: 50-200 µg/ml E. coli: 50-100 µg/ml	3 years at -20°C 2 years at 4°C 3 months at 20-25°C	100 mg/ml	1 g 5 g	ant-hg-1 ant-hg-5
Zeocin™	Cells: 50-300 µg/ml E. coli: 25 µg/ml	2 years at -20°C 1 year at 4°C 6 months at 20-25°C	100 mg/ml	500 mg 1 g 5 g 5 g (bottle) 1 g (powder) 5 g (powder)	ant-zn-05 ant-zn-1 ant-zn-5 ant-zn-5b ant-zn-1p ant-zn-5p
Puromycin	Cells: 1-10 µg/ml E. coli: 100-125 µg/ml	3 years at -20°C 2 years at 4°C 3 months at 20-25°C	10 mg/ml	100 mg 500 mg 500 mg (bottle)	ant-pr-1 ant-pr-5 ant-pr-5b
Phleomycin	Yeast: 10 µg/ml Filamentous Fungi: 25-150 µg/ml	1 year at -20°C 1 year at 4°C 1 month at 20-25°C	20 mg/ml	100 mg 500 mg 250 mg (powder) 500 mg (powder)	ant-ph-1 ant-ph-5 ant-ph-2p ant-ph-5p



Bulk quantity available. For research purposes only. Not intended for use on animals or humans. Zeocin™ is a trademark of InvivoGen.

More information on our website, please visit: www.invivogen.com