

CELL CULTURE

ANTIMICROBIAL AGENTS



Danger is hiding everywhere... keep your cells safe.

- ❖ Potent, fast-acting antimicrobial agents
- ❖ Highly selective: non-toxic to eukaryotic cells
- ❖ Validated through strict quality control

Microbial contamination of cell cultures is a constant and serious threat to your research. Invasive mycoplasmas, bacteria, yeast and fungi can kill or drastically alter your precious cells, leading to erroneous results, lost time and wasted resources.

InvivoGen offers a wide range of highly specific products to help you prevent, detect and eliminate microbial contamination.

Choose from:

Fungin™	Plasmocure™
Normocin™	PlasmoTest™
Normocure™	Primocin™
Plasmocin™	

www.invivogen.com/cell-culture-contamination

Protect your cells

- **Normocin™**

For preventing mycoplasmal, bacterial and fungal infections. A formulation of 3 antibiotics that is routinely added to culture media of animal cells. Normocin™ contains 2 compounds that act on mycoplasmas and on Gram+ and Gram- bacteria by blocking DNA and protein synthesis.

- **Normocure™**

For eliminating Gram+ and Gram- bacterial infections. Unlike common antibiotic treatments such as penicillin/streptomycin (Pen/Strep), Normocure™ is effective against multidrug resistant bacteria. Start to see results in as little as 3 or 4 days, with full elimination by 2 weeks.

- **Primocin™**

For preventing mycoplasmal, bacterial and fungal infections in primary cells. The first formulation to offer complete protection against microbial contaminants. When using Primocin™, there is no need to add penicillin/streptomycin (Pen/Strep).

- **Fungin™**

For preventing and eliminating yeast and fungal infections. A soluble form of Pimaricin that is routinely added to culture media of animal cells. Fungin™ replaces Amphotericin B preparations.

- **PlasmoTest™**

For detecting mycoplasmal infections. Provides a simple, rapid and reliable assay for visual detection of mycoplasmas in cell cultures. PlasmoTest™ is the first assay to utilize cells to signal the presence of mycoplasmas.

- **Plasmocin™ prophylactic™**

For preventing mycoplasmal infections. Contains 2 potent bactericidal components: one that acts on protein synthesis machinery, and one that acts on DNA replication.

- **Plasmocin™ treatment**

For eliminating mycoplasmal infections. Used at a higher concentration than Plasmocin™ prophylactic, it enables full elimination of mycoplasmas in only 2 weeks.

- **Plasmocure™**

For eliminating those rare mycoplasmas that are resistant to Plasmocin™. Combines 2 antibiotics that act through different mechanisms than those in Plasmocin™. A 2-week treatment with Plasmocure™ should be sufficient to completely eliminate the mycoplasmas.

APPLICATION	PRODUCT	MYCOPLASMAS	BACTERIA	YEAST	FUNGI	QANTITY	CAT.CODE
PREVENTION	Plasmocin™ prophylactic	✓	*			25 mg (10 x 1 ml)	ant-mpp
	Normocin™	✓	✓	✓	✓	500 mg (10 x 1 ml)	ant-nr-1
	Primocin™	✓	✓	✓	✓	500 mg (10 x 1 ml)	ant-pm-1
	Fungin™			✓	✓	75 mg (5 x 1.5 ml vials)	ant-fn-1
DETECTION	PlasmoTest™	✓				1 Kit (250 samples)	rep-pt1
ELIMINATION	Plasmocin™ treatment	✓	*			50 mg (2 x 1 ml)	ant-mpt
	Plasmocure™	✓	*			100 mg (1 ml)	ant-pc
	Normocure™		✓			100 mg (2 x 1 ml)	ant-noc
	Fungin™			✓	✓	75 mg (5 x 1.5 ml vials)	ant-fn-1

*Although these products are designed to target mycoplasmas, they are also active against other types of bacteria.

Other pack sizes are available.

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