



**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:

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

[www.invivogen.com/cell-culture](http://www.invivogen.com/cell-culture)

# 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.

## Myc Strip™ NEW

• For immediate detection of mycoplasma contamination in cell cultures. MycoStrip™ is a simple and rapid test based on isothermal PCR. The results are clearly visualized as a band on an immunochromatographic strip.

## PlasmoTest™

• For routine detection of mycoplasma contamination in cell cultures. PlasmoTest™ is a simple and reliable test based on a reporter cellular assay. Colorimetric results are clearly visualized and the amount of contaminants may be quantified using a spectrophotometer.

## Plasmocin™ prophylactic & Plasmocin™ treatment

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

• **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	MYCOPLASMA	BACTERIA	YEAST	FUNGI	QUANTITY	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
	Myc Strip™ NEW	✓				10/20/50/100 tests	rep-mys-x
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](http://www.invivogen.com)