

Fungin™

For the prevention and elimination of fungal contamination in cell cultures

Catalog code: ant-fn-1, ant-fn-2

<http://www.invivogen.com/fungin>

For research use only. Not for human or veterinary use.

Version 18C26-MM

PRODUCT INFORMATION

Content

Fungin™ is supplied as a cell culture tested, sterile filtered light yellow solution at 10 mg/ml. It is available in 2 pack sizes:

- **ant-fn-1:** 5 x 1.5 ml (75 mg)
- **ant-fn-2:** 1 x 20 ml (200 mg)

One 1.5 ml vial is sufficient for 300 ml to 1.5 liters of culture.

One 20 ml bottle is sufficient for 4 and 20 liters of culture.

Shipping and Storage

- Fungin™ is shipped at room temperature. Upon receipt, it can be stored at 4°C or at -20°C. Avoid repeated freeze-thaw cycles.
- The expiry date is specified on the product label.

Note: Product is stable for 2 weeks at room temperature.

QUALITY CONTROL

Each lot is thoroughly tested to ensure the absence of lot-to-lot variation:

- Endotoxin level: < 1 EU/mg
- Physicochemical characterization (pH, appearance)
- Cell culture tested: potency validated on a fungal reference strain

DESCRIPTION

Fungin™ is used as a «routine addition» to eukaryotic cell culture media to prevent fungal (including yeast) contaminations in small or large-scale mammalian cell cultures. Fungin™ can also be used to eliminate fungal (including yeast) contaminations in eukaryotic cell cultures. It is effective against commonly found cell culture contaminants, such as *Candida albicans* and *Aspergillus*.

Fungin™ is an excellent alternative to Amphotericin B, a polyene antifungal antibiotic. Unlike Amphotericin B, Fungin™ is a highly stable compound and it does not need to be dissolved in toxic deoxycholate. Fungin™ provides maximum protection against fungal contamination with minimum cytotoxicity. Fungin™ is cell culture tested, and may be added to media containing commonly used antibacterial agents, such as penicillin and streptomycin (Pen-Strep). Its use has been cited in many publications¹⁻³.

1. Kriek JM. *et al.*, 2016. Female genital tract inflammation, HIV co-infection and persistent mucosal Human Papillomavirus (HPV) infections. *Virology*, 493:247-54.
2. Krutilina R. *et al.*, 2014. MicroRNA-18a inhibits hypoxia-inducible factor 1 α activity and lung metastasis in basal breast cancers. *Breast Cancer Res.*, 16(4):R78.
3. Helias V. *et al.*, 2012. ABCB6 is dispensable for erythropoiesis and specifies the new blood group system Langereis. *Nature Genetics* 44(2):170-3.

COMPOSITION

Fungin™ is a soluble form of Pimaricin, a polyene introduced in the 1950's as an anti-fungal agent. This antimycotic compound kills yeasts, molds and fungi by disrupting ionic exchange through the cell membrane.

METHOD

Preventive use against fungal (including yeast) contamination

To prevent fungal contamination, Fungin™ is used at the concentration of 10 μ g/ml that represents a 1/1000 dilution of the 10 mg/ml stock solution. Refer to the table below to determine the volume of Fungin™ needed to prevent fungal contamination.

Elimination of fungal (including yeast) contamination

Fungin™ treatment requires little hands-on manipulation and is completed in only 5-10 days. It is used at 50 μ g/ml that represents a 1/200 dilution of the 10 mg/ml stock solution. Refer to the table below to determine the volume of Fungin™ needed to eliminate fungal contamination in different volumes of growth medium.

Fungin™ use	T25 with 5 ml medium	T75 with 15 ml medium	500 ml bottle
For prevention	5 μ l	15 μ l	500 μ l
For decontamination	25 μ l	75 μ l	2.5 ml

RELATED PRODUCTS

Product	Description	Cat. Code
Normocin™	Antimicrobial agent	ant-nr-1
Normocure™	Antibacterial agent	ant-noc
Plasmocin™ Prophylactic	Anti-mycoplasma agent	ant-mpp
Plasmocin™ Treatment	Mycoplasma removal agent	ant-mpt
PlasmoTest™	Mycoplasma detection kit	rep-pt1
Primocin™	Antimicrobial for primary cells	ant-pm-1

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: info@invivogen.com