

## Instruction Manual

Product Name	Product Description	Size	Catalog Number
BIOMYC-2 Antibiotic Solution, 100X	BIOMYC-2 Antibiotic Solution, 100X	10 ml	PK-CC03-037-1D
		20 ml	PK-CC03-037-1C
		100 ml	PK-CC03-037-1B

### Introduction

The contamination of cells with mycoplasma is a very common problem, even though it often remains unnoticed since no cloudiness appears in the cell culture. Nevertheless, the contamination often causes biochemical changes as well as changes in the immunological properties of the cells. Since mycoplasma-infected cells cannot always be discarded, many complicated methods have been suggested for the elimination of the mycoplasma.

PromoKine is now offering a combination of antibiotics, which have been shown to be effective in the elimination of mycoplasma species that account for 90% of the contamination found in cell cultures. When used according to the following instructions, no cytotoxic effects will occur.

### Product Description

BIOMYC-1 is based on the antibiotic tiamutin, which is produced by the fungus *pleurotus mutilus*. BIOMYC-2 is based on minocycline, which is a tetracycline derivative. These two antibiotic solutions are generally used sequentially in combination.

### Storage & Stability

Store at -20°C.

### Instructions for Use

1. Do not use the two solutions together, but rather sequentially.
2. Add 1 ml BIOMYC-1 to 100 ml medium, and keep the contaminated cells in this mixture for 4 days. Any fresh medium added should also contain BIOMYC-1.
3. After 4 days, add 1 ml BIOMYC-2 to 100 ml fresh medium, and maintain the cells in this second mixture for 3 days.
4. The above, together, are considered as one treatment cycle. It may be necessary to repeat this cycle 2-3 times.
5. During the process, the cells can be tested for mycoplasma contamination, and results can then be used to shorten the process when possible.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.