

Instruction Manual

Product Name	Product Description	Size	Catalog Number
BIOMYC-3 Antibiotic Solution, 100X	BIOMYC-3 Antibiotic Solution, 100X	10 ml	PK-CC03-038-1D
		20 ml	PK-CC03-038-1C
		100 ml	PK-CC03-038-1B

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-3 is based on the ciprofloxacin antibiotic, which is a member of the fluoroquinolone group. Many mycoplasma species have been found to be sensitive to BIOMYC-3, including *A.laidlawii*, *M. orale*, *M. hyorhinae*, *M. fermentans*, and *M. arginini*. These species are responsible for most of the contamination in cell culture. At the concentrations recommended for use, no cytotoxic effects have been found, and the treatment is quite easy to perform.

Storage & Stability

Store at -20°C.

Instructions for Use

1. Add 1 ml BIOMYC-3 to 100 ml medium.
2. Continue the treatment for a total of 14 days, while changing the medium (containing BIOMYC-3) every 2-3 days.
3. Retain the cells in the growth medium for an additional 14 days before re-testing for mycoplasma.

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