

Mycoplasma Cleaner

Catalogue number	Presentation
MB46401	250 mL
MB46402	500 mL (Refill)

Features

- Ready-to-use solution specifically designed to decontaminate laboratory surfaces and apparatus.
- It offers a wide spectrum of activity, effectively targeting *Mycoplasma spp.*, *Staphylococcus aureus*, *Streptococcus faecalis*, *Escherichia coli*, *Proteus mirabilis*, *Pseudomonas aeruginosa*, *Candida albicans*, as well as HBV/HIV, *Papovavirus*, *Adenovirus*, *Poliovirus*, and other contaminant microorganisms.
- The cleaner is conveniently packaged in a user-friendly, lightweight bottle equipped with a trigger spray for easy application.
- Using this cleaner simplifies the decontamination process, saving both time and effort.
- A refill option is available for added convenience.

Description

Mycoplasma Cleaner is a highly effective decontaminating solution with broad-spectrum activity, designed to target not only Mycoplasma but also viruses, bacteria, and fungi. This solution provides a safe and reliable approach to decontamination in laboratory settings, particularly on benches, safety cabinets, incubators, cell storage boxes, and liquid nitrogen containers, among other surfaces and equipment.

Mycoplasma species, which are common contaminants in cell cultures, can be challenging to detect using standard microscopy techniques due to their small size. As a result, Mycoplasma infections often go unnoticed, leading to cell infection and a range of detrimental effects. Contamination can occur through media spills from contaminated cell cultures during loading and unloading procedures in biosafety cabinets, resulting in surface contamination where Mycoplasma can survive for extended periods. To prevent cross-contamination, it is crucial to promptly isolate and eliminate infected cultures, as even low levels of Mycoplasma infection can disrupt cell growth, alter metabolism, and protein synthesis. Regular and thorough cleaning and disinfection of all surfaces and equipment are vital measures to combat and prevent Mycoplasma contamination in the laboratory.

By utilizing the Mycoplasma Cleaner and implementing proper hygiene practices, you can effectively mitigate the risk of mycoplasma contamination, ensuring the integrity of your cell cultures and the reliability of your research.

As an additional measure for efficient contamination control, we highly recommend conducting routine testing of cell culture materials using our **Mycoplasma spp. qPCR Kit (Cat. No. MD05511)**. This kit is specifically designed for the in vitro detection of *Mycoplasma spp.* genomes, providing a fast, simple, and effective method to identify mycoplasma-contaminated cell cultures. The real-time qPCR-based method utilized by the *Mycoplasma spp.* qPCR Kit can be conveniently performed in any standard molecular biology laboratory, offering a valuable tool to ensure the integrity and reliability of your cell culture experiments.

The Mycoplasma Cleaner has undergone extensive testing and validation to ensure optimal performance in diverse laboratory settings. Its effectiveness is scientifically proven and supported by robust evidence, providing you with the assurance you need for critical decontamination requirements.

Shipping & Storage Conditions

This product is shipped at room temperature. Upon receipt, store it at 15 °C to 25 °C.

Standard Protocol

Use gloves as personal protective equipment when manipulating this product.

Decontamination of surfaces:

1. Spray Mycoplasma Cleaner directly onto the contaminated surfaces until they are completely moistened.
2. Allow the product to act for 5 minutes.
3. Gently wipe the surface with a paper towel until dry. The Cleaner does not leave any residue.

Note: Prior to use, it is recommended to perform a surface compatibility test as Mycoplasma Cleaner may cause discoloration on sensitive surfaces.

Decontamination of apparatus:

1. Apply Mycoplasma Cleaner onto a paper towel.
2. Vigorously rub all surfaces of the apparatus, ensuring thorough coverage.
3. Thoroughly dry the surface with a clean paper towel.

Please note that while our recommendations provide a robust decontamination solution, we encourage users to verify the suitability of Mycoplasma Cleaner for their specific purposes and applications. Do not add the cleaner to the reaction.

FOR REFERENCE USE ONLY

For life science research only. Not for use in diagnostic procedures.