

April 2020

Keywords or phrases:

Mycoplasma; pipetting; decontamination; cell culture; filter tips; avoiding contamination

Minimize the risk of mycoplasma contamination in your cell cultures

Understand the source and prevention

Emilia Varhimo¹, Sandra Söderholm²

1. Pipette and Dispensing, Sartorius Biohit Liquid Handling Oy, Helsinki, Finland

2. Lab Essentials Applications Development, Sartorius Biohit Liquid Handling Oy, Helsinki, Finland

Correspondence:

E-Mail: lhinfo.Finland1@Sartorius.com

Abstract

Preventing mycoplasma contamination in the cell culture lab can be a challenge. However, you can significantly reduce the risk of mycoplasma contamination and spread in your cell culture lab with good cell culture practices, by selecting easy-to-clean pipettes and sterility promoting consumables, and by regularly testing.

Find out more: www.sartorius.com/en/products/pipetting-and-dispensing

Introduction

Mycoplasma (shown in Figure 1) contamination is prevalent in cell culture labs: in one study, researchers identified mycoplasma contamination in 11% of the 10,000 cell lines they tested (Olarerin-George et al. 2015). Part of the reason mycoplasma contamination is so prevalent is that mycoplasma spreads efficiently even in routine subculturing. In another study, researchers detected live mycoplasma outside of the flask, on a hemocytometer, on the pipette, and outside of the pipette discard pan after working with mycoplasma-infected cell culture in a laminar flow hood. Indeed, live mycoplasma were recovered from the surface of the laminar flow hood even four to six days later. And a clean culture that was subcultured once a week in the same hood following work with the contaminated cells tested positive for mycoplasma after only 6 weeks (reviewed in Nikfarjam et al 2012, original article McGarrity GJ (1976).

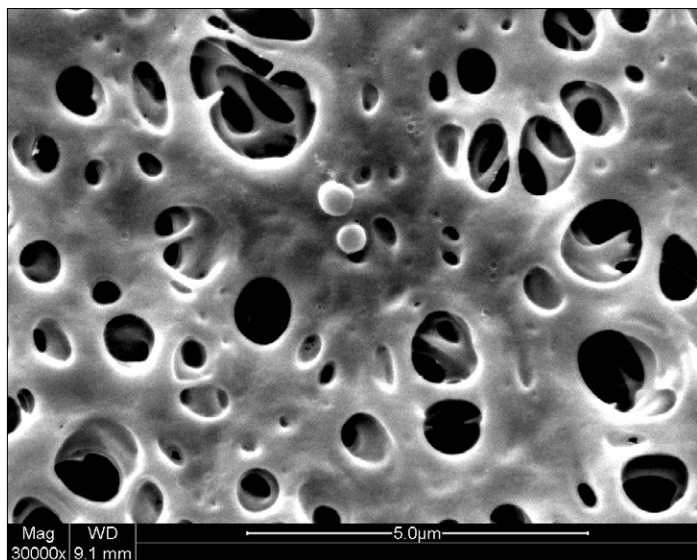


Figure 1. Scanning electron microscope image of mycoplasma (*A. laidlawii*) retained on 0.1 μ m PES membrane.

What are mycoplasma?

Mycoplasma are a genus of bacteria that lack a cell wall. Because they lack a cell wall, they can take on different shapes, from round to elongated, and they are resistant to antibiotics (Bebear et al. 2011). Due to their small size (0.2 – 0.8 μ m), mycoplasma can penetrate sterilizing-grade filters, making them a very difficult contaminant to eliminate. Their small size also makes them difficult to spot within cell cultures (Masters et al. 2007; Volokhov et al. 2008).

In cell culture, mycoplasma inhibit protein biosynthesis and cell growth and alter RNA and DNA synthesis, thus they can affect research results. In cell based therapeutics and in advanced therapy medical products (ATMPs), mycoplasma contamination can cause immune reactions, chromosomal aberrations, or altered proliferation characteristics, making therapy unsafe for use in patients (Martins et al. 2014).

Potential routes for mycoplasma contamination include:

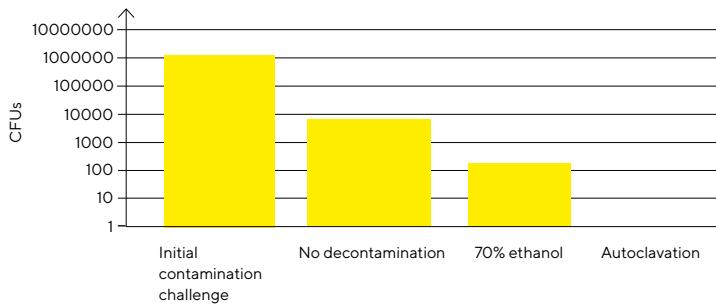
- Contaminated media, reagents or instruments
- Laboratory personnel (approximately 80.6% of technicians are mycoplasma carriers)
- Infected cultures obtained from other labs

Mycoplasma persist on pipette surfaces

Pipettes are frequently exposed to contaminants and potentially can carry the contamination forward. For this reason, most pipettes are made to be either fully autoclavable or have dispensing head parts that can be autoclaved. We performed a study to compare the efficiency of decontamination by autoclaving and wiping with ethanol. We contaminated the surface of three pipettes with 10^6 mycoplasma CFUs on defined areas and then left them in a laminar flow for 24 hours. After this the pipettes were either autoclaved or wiped with ethanol. We recovered no mycoplasma after autoclaving, compared to about 100 CFU after wiping with 70% ethanol (see Figure 2). Without any decontamination procedure, about 10 000 CFUs could be recovered from pipette after 24 hours, indicating the persistence of mycoplasma.

This result recapitulates the results in a similar study by Eterpi et al (2010). They found that decontamination with 70% ethanol reduced the number of recovered bacteria by more than one log. They also found that autoclaving pipettes completely removes mycoplasma contamination.

To mitigate the risk of mycoplasma contamination, use autoclavable pipettes for cell culture work. Sartorius's fully autoclavable Tacta pipette is a good choice for all cell culture work.



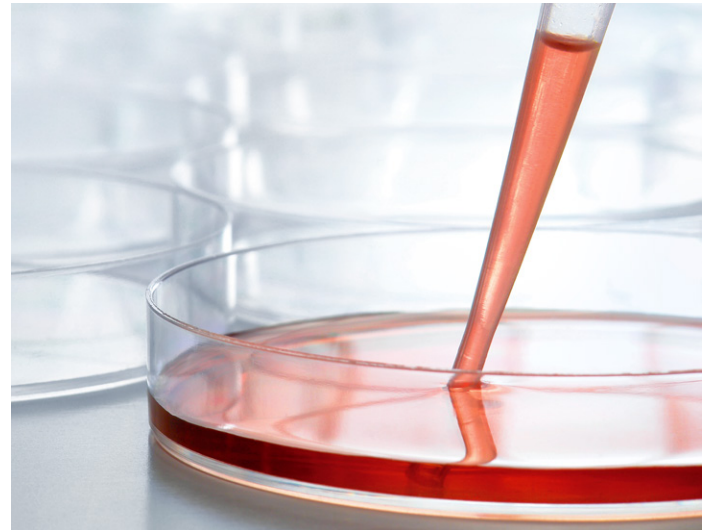
Best practices to avoid mycoplasma contamination

You can achieve contamination-free cell cultivation with these practical procedures:

1. Implement Good Cell Culture Practice

Good Cell Culture Practice (Coecke S, 2005) is a set of general cell culture guidelines to prevent contamination without needing to resort to the use of antibiotics. Routine use of antibiotics in cell culture media is not recommended. Not only are antibiotics often ineffective against mycoplasma, but they also promote antibiotic resistance and can hide low-level persistent contamination (Drexler & Uphoff, 2002). In addition, antibiotics in media may have an effect on the outcome of cell culture experiments.

Key guidelines in Good Cell Culture Practice include: cultivation of only one cell line at a time; clearly marking the pipettes, pipetting controllers, and pipette tips used for cell culture work; only using these pipettes and tips in the cell culture laboratory; and never moving them from the cell culture laboratory to another and back again. Only use reagents designated exclusively for cell culture work, and do not use stock solutions designated for other applications.



2. Select easy-to-clean and autoclavable pipettes for cell culture work

Maintenance of pipettes and liquid handling controllers play an important part in reducing risks of spilling and splashing of liquids, a key source of contamination. Laboratories should set guidelines and protocols for regular decontamination and maintenance. According to Eterpi et al. (2010), effective cleaning methods for removing bacterial contaminations are:

- Autoclaving
- Alkaline cleaners
- Ethanol Vaporized hydrogen peroxide VHP

There are significant differences between pipette designs that affect their cleaning protocols. For instance, depending on the design, a disassembled pipette could be between three or more than twenty parts. To make cleaning and sterilizing easy and reliable, choose pipetting tools that are fully autoclavable, easy-to-clean, and tolerate effective cleaning methods, such as the Sartorius Tacta pipette.





3. Select sterile filter tips with protective wrappings

Filter tips are the safest choice for cell culture work, as they provide the best protection for both the samples and the pipette and they prevent mycoplasma spreading by aerosols. Autoclaving is not suitable for filter tips because the filter material, polyethylene, does not tolerate high temperatures. Therefore, pre-sterilized pipette tips are most commonly sterilized with electron beam (beta) irradiation. Sartorius uses e-beam pre-sterilization with a sterility assurance level (SAL) of 10^{-6} . This means that there is a maximum probability of one viable microorganism in 1 million sterilized items. This SAL level is accepted for pharmacopoeial sterilization procedures. Sartorius evaluates the e-beam radiation dose four times a year with independent dose audits.

Dust and dirt collects on the surface of tip boxes when storing tip racks in warehouses and lab shelves. Therefore, choose individually wrapped tip boxes, because the wrapping ensures sterility during storage. Only open the wrapping just before use and put the unwrapped box straight into the laminar flow cabinet to minimize carryover of dirt to the clean working area.



4. Test regularly

Test all cell lines regularly for mycoplasma contamination, especially new cell lines. Keep new cell lines isolated in a separate incubator until they can be reliably shown free of mycoplasma. See Table 1 for some of the common mycoplasma detection methods, and their advantages and disadvantages. Many researchers prefer using mycoplasma PCR/qPCR kits because they are very sensitive and results can be obtained fast. As PCR/qPCR cyclers are common in most laboratories, the testing is easy to conduct.



Detection limits in different cell types

Method	Advantages	Disadvantages
Microbial culture	+ High sensitivity and specificity	<ul style="list-style-type: none"> Requires special microbiology lab Slow (takes up to 28 days)
Direct DNA stain (DAPI, Hoechst)	+ Rapid and cheap	<ul style="list-style-type: none"> Interpretation of results is difficult and subjective Sensitivity is very low
Mycoplasma kit (PCR/qPCR)	+ Rapid and very sensitive + Easy to interpret results	<ul style="list-style-type: none"> No discrimination between live and dead bacteria

Table 1. Mycoplasma testing methods

5. Filter with 0.1 µm filter

Autoclaving is an efficient way to kill mycoplasma, but unsuitable for some media and reagents, because heat destroys many nutrients and growth factors. Standard sterile filtration with pore size of 0.22-µm that is generally efficient in preventing the passage of bacteria and fungi is inadequate for mycoplasma that have a cell size of 0.2-0.8 µm and can pass through pores of > 0.1 µm in size. Therefore, to prevent mycoplasma contamination, use the 0.1 µm pore size instead of standard 0.22 µm sterile filters when filtering cell culture reagents and media (Roche et al. 1992).



References

Bebear C, Pereyre S, Peuchant O. Mycoplasma pneumoniae: susceptibility and resistance to antibiotics. *Future Microbiol* (2011); 6; 423-431.

Coecke S, Balls M, Bowe G, Davis J, Gstraunthaler G, Hartung T, Hay R, Merten O-W. Guidance on Good Cell Culture Practice. *ATLA* (2005); 33; 261-287.

Drexler HG & Uphoff CC. Mycoplasma contamination of cell cultures: Incidence, sources, effects, detection, elimination, prevention. *Cytotechnology* (2002); 39; 75-90.

Eterpi M, McDonnell G, Thomas V. Decontamination efficacy against Mycoplasma. *Letters in Applied Microbiology* (2010); 52; 150-155.

Martins JP, Santos JM, de Almeida JM, Filipe MA, de Almeida MV, Almeida SC, Água-Doce A, Varela A (2014). Towards an advanced therapy medicinal product based on mesenchymal stromal cells isolated from the umbilical cord tissue: quality and safety data. *Stem Cell Res Ther* (2014); 5(1); 9.

Masters JR, Stacey GN. Changing medium and passaging cell lines. *Nat. Protoc.* (2007); 2; 2276-2284.

McGarrity GJ. Spread and control of mycoplasma infection of cell cultures. *In vitro* (1976); 12 (9); 643-648.

Nikfarjam L & Farzaneh P. Prevention and Detection of Mycoplasma Contamination in Cell Culture. *Cell J* (2012); 13(4); 203-212.

Olarerin-George AO & Hogenesch JB. Assessing the prevalence of mycoplasma contamination in cell culture via a survey of NCBI's RNA-seq archive. *Nucl Acid Res* (2015); 43(5); 2535-2542.

Roche KL, Levy RV. Methods used to validate microporous membranes for the removal of Mycoplasma. *Biopharm* 1992;22-33.

Volokhov DV, Kong H, George J, Anderson C, Chizhikov VE. Biological enrichment of mycoplasma agents by cocultivation with permissive cell cultures. *Appl. Environ. Microbiol.* (2008); 74; 5383-5391.

Sales and Service Contacts

For further contacts, visit
www.sartorius.com

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

Finland

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki
Finland
Email: lhinfo.Finland1@Sartorius.com