

# SARTORIUS



## Rapid Mycoplasma Detection

Microsart® AMP Extraction

Microsart® ATMP Mycoplasma

A cutting-edge solution for the reliable and efficient detection of Mycoplasma contaminations

- Microsart® AMP Extraction for efficient extraction of Mycoplasma DNA
- Microsart® ATMP Mycoplasma for real-time PCR

# Rapid Mycoplasma Detection

Mycoplasma contamination in cell cultures is a serious concern in many applications, as it can lead to unreliable results and wasted time and resources. To combat this quality issue, Microsart® ATMP Mycoplasma offers a reliable and easy-to-use solution for detecting mycoplasma contaminations in cell culture samples.

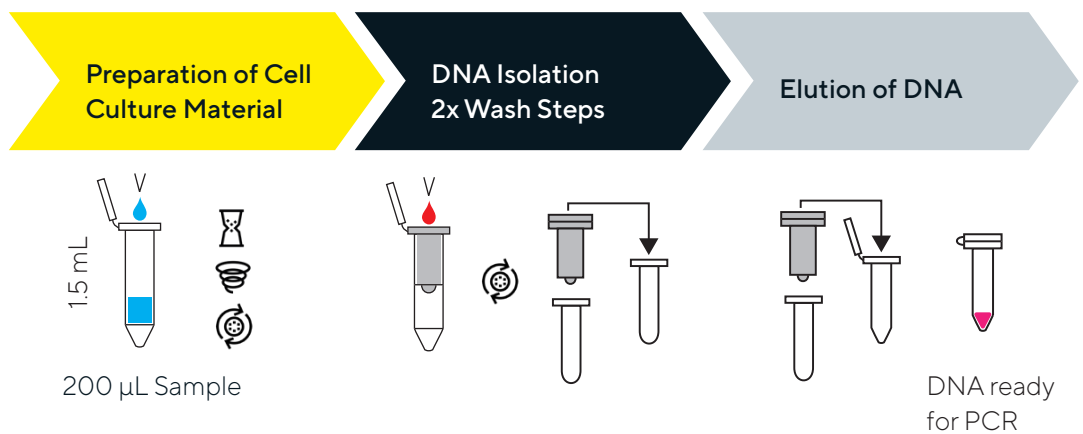
Based on real-time PCR technology Microsart® ATMP Mycoplasma allows for highly sensitive and specific detection of mycoplasma DNA. It is validated to detect Mycoplasma contaminations of  $\leq 10$  CFU/mL.

One of the key advantages of the Microsart® ATMP Mycoplasma kit is its rapid time-to-result and its simplicity of use. Using our Microsart® AMP Extraction and Microsart® ATMP Mycoplasma kits is a simple two-step process: DNA isolation and amplification. The kit includes all necessary reagents and controls to ensure accurate and reliable results, ruling out false positives and negatives. The extraction process is especially adapted for efficient isolation of Mycoplasma DNA.

The Microsart® ATMP Mycoplasma kits offer a reliable, user-friendly solution for the rapid detection of mycoplasma contaminations in cell culture samples. With its proven performance and ease of use, it is an essential tool for USP/EP conform mycoplasma release testing.

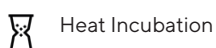
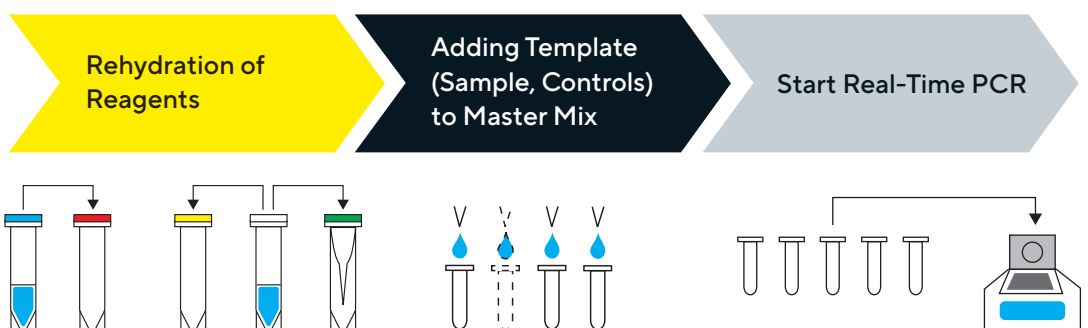
## DNA Extraction

A column-based extraction designed to capture all DNA of fragile Mycoplasma



## PCR Setup

Uniform in all Microsart® ATMP kits



# Microsart® real-time PCR Kits and PCR related products

## Bacteria & Fungi Detection

Order No.	Description	Quantity
SMB95-1007	Microsart® ATMP Sterile Release	10 patient sample releases
SMB95-1008	Microsart® ATMP Bacteria	100 reactions
SMB95-1012	Microsart® ATMP Fungi	100 reactions
SMB95-1009	Microsart® Research Bacteria	25 reactions
SMB95-1014	Microsart® Research Fungi	25 reactions

## DNA Extraction Kit

Order No.	Description	Quantity
SMB95-2001	Microsart® ATMP Extraction	Reagents for 50 extractions
SMB95-2003	Microsart® AMP Extraction (only for Mycoplasma qPCR)	Reagents for 50 extractions

## Mycoplasma Detection Kits for qPCR

Order No.	Description	Quantity
SMB95-1001 1002	Microsart® AMP Mycoplasma	25   100 reactions
SMB95-1003 1004	Microsart® ATMP Mycoplasma	25   100 reactions
SMB95-1005	Microsart® Research Mycoplasma	25 reactions

## Microsart® Validation Standard according to EP 2.6.7 and USP <63> for Mycoplasma species and EP 2.6.1, EP 2.6.27 and USP <71> for other bacteria and fungi

3 vials with 10 CFU/vial for Mycoplasma species and 6 vials with 99 CFU/vial for other bacteria and all fungi

Order No.	Description
SMB95-2005	<i>Bacillus subtilis</i>
SMB95-2006	<i>Pseudomonas aeruginosa</i>
SMB95-2007	<i>Kocuria rhizophila</i>
SMB95-2008	<i>Clostridium sporogenes</i>
SMB95-2009	<i>Bacteroides vulgatus</i>
SMB95-2010	<i>Staphylococcus aureus</i>
SMB95-2011	<i>Mycoplasma arginini</i>
SMB95-2012	<i>Mycoplasma orale</i>
SMB95-2013	<i>Mycoplasma gallisepticum</i>
SMB95-2014	<i>Mycoplasma pneumoniae</i>
SMB95-2015	<i>Mycoplasma synoviae</i>
SMB95-2016	<i>Mycoplasma fermentans</i>
SMB95-2017	<i>Mycoplasma hyorhinis</i>
SMB95-2018	<i>Acholeplasma laidlawii</i>
SMB95-2019	<i>Spiroplasma citri</i>
SMB95-2020	<i>Mycoplasma salivarium</i>

Order No.	Description
SMB95-2037	<i>Candida albicans</i>
SMB95-2038	<i>Aspergillus brasiliensis</i>
SMB95-2039	<i>Aspergillus fumigatus</i>
SMB95-2040	<i>Penicillium chrysogenum</i>
SMB95-2041	<i>Candida glabrata</i>
SMB95-2042	<i>Candida krusei</i>
SMB95-2043	<i>Candida tropicalis</i>

## Microsart® Calibration Reagent

1 vial, 10<sup>8</sup> genomes/vial for all bacteria and 10<sup>6</sup> genomes/vial for all fungi


Order No.	Description
SMB95-2021	<i>Mycoplasma arginini</i>
SMB95-2022	<i>Mycoplasma orale</i>
SMB95-2023	<i>Mycoplasma gallisepticum</i>
SMB95-2024	<i>Mycoplasma pneumoniae</i>
SMB95-2025	<i>Mycoplasma synoviae</i>
SMB95-2026	<i>Mycoplasma fermentans</i>
SMB95-2027	<i>Mycoplasma hyorhinis</i>
SMB95-2028	<i>Acholeplasma laidlawii</i>
SMB95-2029	<i>Spiroplasma citri</i>
SMB95-2030	<i>Bacillus subtilis</i>
SMB95-2031	<i>Pseudomonas aeruginosa</i>
SMB95-2032	<i>Kocuria rhizophila</i>
SMB95-2033	<i>Clostridium sporogenes</i>
SMB95-2034	<i>Bacteroides vulgatus</i>
SMB95-2035	<i>Staphylococcus aureus</i>
SMB95-2036	<i>Mycoplasma salivarium</i>
SMB95-2044	<i>Candida albicans</i>
SMB95-2045	<i>Aspergillus brasiliensis</i>
SMB95-2046	<i>Aspergillus fumigatus</i>
SMB95-2047	<i>Penicillium chrysogenum</i>
SMB95-2048	<i>Candida glabrata</i>
SMB95-2049	<i>Candida krusei</i>
SMB95-2050	<i>Candida tropicalis</i>

## 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

 For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)