

SARTORIUS



Rapid Solutions

Microsart® ATMP Bacteria, Fungi and Mycoplasma

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

- Microsart® ATMP Mycoplasma
- Microsart® ATMP Bacteria & Microsart® ATMP Fungi PCR
- Microsart® ATMP Sterile Release for single sample release

Rapid Solutions

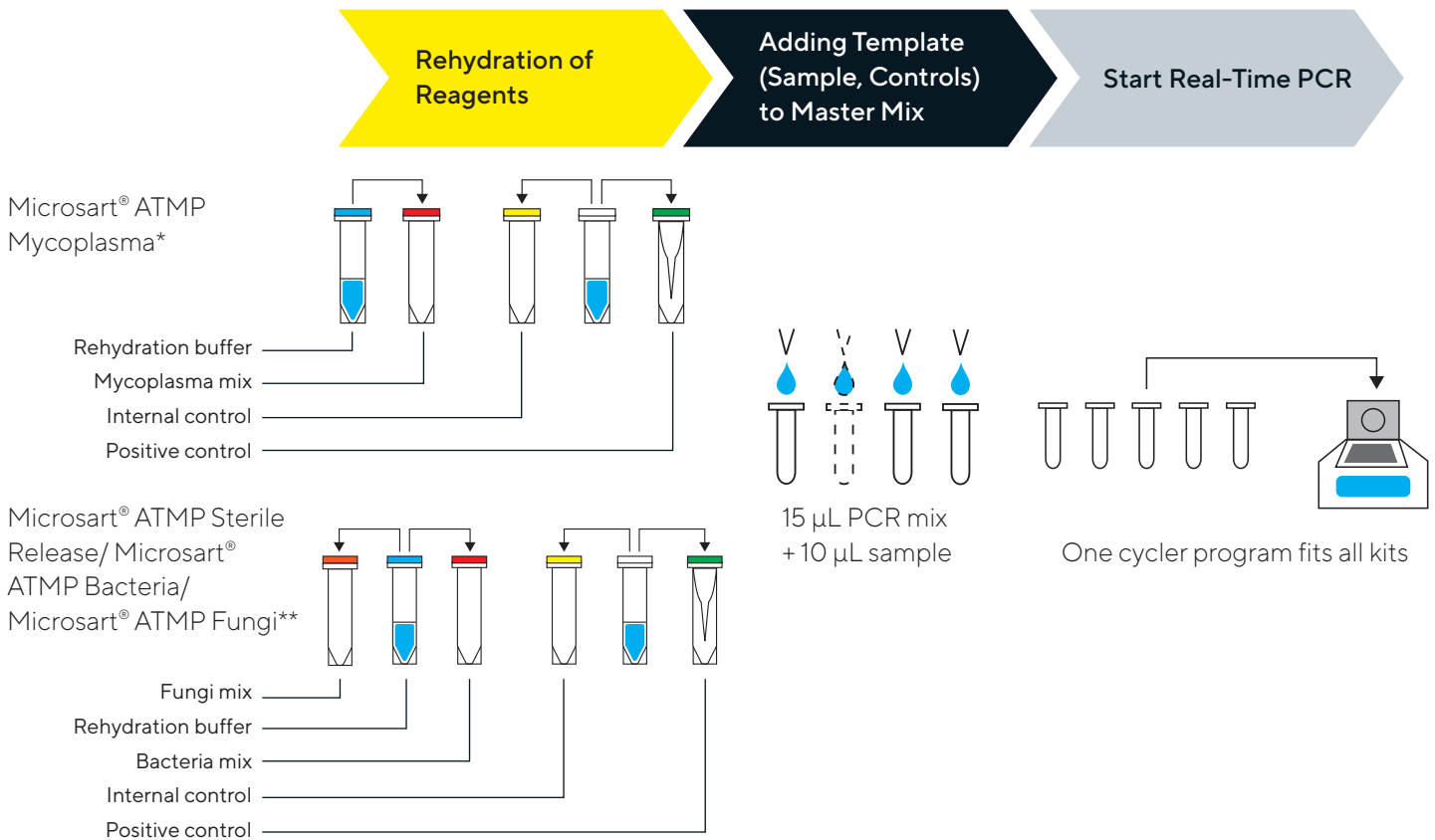
Sartorius Microsart ATMP® Mycoplasma, Microsart® ATMP Bacteria, and Microsart® ATMP Fungi are essential tools for the pharmaceutical industry to ensure the quality of Advanced Therapy Medicinal Products (ATMPs) or other short-live products during production.

Mycoplasma, bacteria, and fungi contamination in ATMPs can have serious consequences for patient safety and process efficiencies. Microsart® ATMP Mycoplasma detects Mycoplasma DNA using a highly specific and sensitive real-time PCR technology, while Microsart® ATMP Bacteria and Microsart® ATMP Fungi detect bacterial and fungal DNA, respectively.

One of the key advantages of these kits is their time-to-result, sensitivity and specificity. All three kits share a common protocol and can be used in line to simultaneously detect Mycoplasma, bacteria & fungi in one sample. 15 µL of each master mix is mixed with 10 µL of the sample DNA, positive control DNA or negative control. This streamlined process simplifies laboratory workflows and reduces the potential for errors, making the kits a popular choice for quality control laboratories in the pharmaceutical industry.

Microsart® ATMP Mycoplasma, Microsart® ATMP Bacteria, and Microsart® ATMP Fungi are essential tools for quality control in pharmaceutical production, enabling pharmaceutical manufacturers to meet regulatory requirements and ensure the safety and efficacy of ATMPs for patients.

PCR Setup



* The Microsart® ATMP Mycoplasma Kit is validated with the Microsart® AMP Extraction kit

** The Microsart® ATMP Bacteria/Fungi kits are validated with the Microsart® ATMP Extraction kit

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, 10⁸ genomes/vial for all bacteria and 10⁶ 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