

Microsart® Validation Standard

Prod. No. SMB95-2011	Mycoplasma arginini
Prod. No. SMB95-2012	Mycoplasma orale
Prod. No. SMB95-2013	Mycoplasma gallisepticum
Prod. No. SMB95-2014	Mycoplasma pneumoniae
Prod. No. SMB95-2015	Mycoplasma synoviae
Prod. No. SMB95-2016	Mycoplasma fermentans
Prod. No. SMB95-2017	Mycoplasma hyorhinis
Prod. No. SMB95-2018	Acholeplasma laidlawii
Prod. No. SMB95-2019	Spiroplasma citri
Prod. No. SMB95-2020	Mycoplasma salivarium

10 CFU

For use in research and quality control

Manufactured by:



Symbols

LOT

Lot No.

REF

Order No.



Expiry date



Store at



Content

Contents

1. Intended Use	5
2. Explanation of the Product	5
3. Notes on the Test Procedure	6
4. Reagents	7
5. Needed but not included	8
6. Precautions	9
7. Test Procedure	9
8. Related Products	10

1. Intended Use

Microsart® Validation Standard is used to validate the robustness and sensitivity of NAT-based mycoplasma detection methods in combination with cell cultures, cell culture media components, and cell culture-derived biologicals (e.g. Advanced Therapy Medicinal Products, ATMPs, such as cell autologous transplants), according to the European Pharmacopoeia 2.6.7 (EP 2.6.7 „Mycoplasmas“).

2. Explanation of the Product

In a regulated environment, microbial detection can be extremely time-consuming if it relies on conventional culture methods. Rapid microbial tests such as NAT-based methods (e.g. PCR) are formally valid alternatives to traditional culture methods, only upon a comprehensive assay validation. In this regard, the EP 2.6.7 requires the validation of the test sensitivity and robustness with respect to the sample matrix and lab precision. In addition, the applied analytical method shall show a performance equal or better than the compendial culture procedure. As culturing living mycoplasma for validation purposes represents a highly demanding when not impracticable task for most laboratories, safe and easy-to-use alternatives are needed.

Microsart® Validation Standards are inactivated and therefore non-infectious mycoplasma preparations. They are titrated to 10 Colony-forming Units/ml (CFU/ml), the sensitivity limit indicated for NAT-based methods like PCR to replace the traditional culture method.

The following species are available:

Mycoplasma arginini	ATCC 23838; NCTC 10129
Mycoplasma orale	ATCC 23714; NCTC 10112; DSM 25590
Mycoplasma gallisepticum	ATCC 19610; NCTC 10115; DSM 19817
Mycoplasma pneumoniae	ATCC 15531; NCTC 10119; DSM 22911
Mycoplasma synoviae	ATCC 25204; NCTC 10124; DSM 21430
Mycoplasma fermentans	ATCC 19989; NCTC 10117
Mycoplasma hyorhinis	ATCC 17981; NCTC 10130; DSM 25591
Acholeplasma laidlawii	ATCC 23206; NCTC 10116; DSM 23060
Spiroplasma citri	ATCC 27556; NCTC 10164; DSM 21846
Mycoplasma salivarium	ATCC 23064; NCTC 10113

The mycoplasma used for the manufacture of Microsart® Validation Standard are low passage reference strains cultivated in culture broth (Hayflick and Frey medium), as described in EP 2.6.7. The cultures are harvested in the logarithmic growth phase to avoid a high ratio of dead mycoplasma particles, titrated in culture broth and plated on Hayflick and Frey medium for quantification based on CFU.

Each vial contains inactivated mycoplasma particles corresponding to 10 CFU. The relevant sample matrix can be added directly into the vial. The derived sample is expected to be tested positive by a valid NAT-based assay. The inactivated mycoplasma preparation cannot be used for the culture method. For maximal sensitivity, the mycoplasma DNA should be extracted prior to PCR. For this purpose, we developed Microsart® AMP Extraction (Sartorius Prod. No. SMB95-2003). For PCR, we recommend using Microsart® AMP Mycoplasma (SMB95-1001/ SMB95-1002), Microsart® ATMP Mycoplasma (Sartorius Prod. No. SMB95-1003/ SMB95-1004), or Microsart® Research Mycoplasma (Sartorius Prod. No. SMB95-1005/1006).

3. Notes on the Test Procedure

1. This leaflet must be fully understood in order to successfully use Microsart® Validation Standard. The reagents should not be used beyond their shelf life.
2. Any deviation from the described method can affect the results.
3. Inhibition of PCR may be caused by the sample matrix. Negative controls should be processed with the same sample matrix.
4. For each sample matrix, at least one negative control should be tested. Resulting Ct values can be compared to lot-specific Ct values specified in the respective Certificate of Analysis.
5. Participation in external quality control programs, such as those offered by Minerva Biolabs GmbH (www.minerva-biolabs.com), is recommended.

4. Reagents

Each product contains 3 vials of mycoplasma particles as well as 2 vials containing the same carrier matrix as the mycoplasma vials for the preparation of corresponding negative controls. All components are lyophilized for maximal product stability.

All particles have been inactivated prior to lyophilization. The expiry date of the unopened package is specified on the package label. The kit components are stored at +2 to +8 °C upon arrival, at low humidity conditions.

The lot specific Certificates of Analysis can be downloaded from the MySartorius portal (<https://my.sartorius.com>).

Component Label Information	Order No.	Quantity	Cap Color
Mycoplasma			
Acholeplasma	SMB95-2011-	3 × lyophilized	green
Spiroplasma	SMB95-2020		
Negative Control		2 × lyophilized	white

5. Needed but not included

Microsart® Validation Standard contains the positive and negative preparations to perform the test. General industrial supplies and reagents, usually available in PCR laboratories are not included:

Consumables

- Laboratory gloves
- PCR reaction tubes for the specific qPCR device
- PCR grade pipette filter tips

Equipment

- PCR cycler
- Microcentrifuge for 1.5 ml reaction tubes
- Vortex
- Rack for 1.5 ml tubes
- Pipettes (Sartorius)
 - mechanical
 - 0.5 – 10 µl Sartorius Prod. No. LH-729020
 - 10 – 100 µl Sartorius Prod. No. LH-729050
 - 100 – 1000 µl Sartorius Prod. No. LH-729070
 - or electrical
 - 0.2 – 10 µl Sartorius Prod. No. 735021
 - 10 – 300 µl Sartorius Prod. No. 735061
 - 50 – 1000 µl Sartorius Prod. No. 735081

For DNA extraction and PCR analysis, the following kits are required additionally:

- Mycoplasma DNA extraction system. We recommend the Microsart® AMP Extraction (Prod. No. SMB95-2003).
- Mycoplasma DNA PCR detection system. We recommend the Microsart® AMP Mycoplasma (Prod. No. SMB95-1001/1002), Microsart ATMP Mycoplasma (Prod. No. SMB95-1003/1004) or Microsart® Research Mycoplasma (Prod. No. SMB95-1005/1006).

6. Precautions

For *in vitro* use in research and quality control. This product should be used only by trained persons. All samples should be considered potentially infectious and handled according to the local or national regulations. This product does not contain hazardous substances and may be disposed of according to local guidelines.

7. Test Procedure

-
1. Centrifuge the tube(s) briefly to collect the lyophilized material at the bottom.

 2. Add 1 ml of the sample matrix of interest to each vial.

 3. Incubate for at least 5 minutes at room temperature. Proceed immediately when the lyophilized CFU material is completely dissolved.

 4. Vortex for 10 sec and optionally spin down 5 sec to remove droplets from the lid.

 5. Use the volume of sample required by the selected sample preparation kit. After DNA extraction, proceed to PCR.

All reagents and samples must be equilibrated to room temperature before use.

It is highly recommended to perform suitable DNA extraction of the samples prior to PCR in order to reduce the risk of PCR inhibition and maximize sensitivity.

The Negative Control vials contain exactly the same components (carrier matrix) as the Mycoplasma vials except for the mycoplasma particles. For a valid interpretation of the test results, the Negative Controls should be rehydrated with the sample matrix of interest and processed in parallel to the samples, in a suitable number of replicates.

8. Related Products

Detection Kits for qPCR

SMB95-1001/1002	Microsart® AMP Mycoplasma	25/100 tests
SMB95-1003/1004	Microsart® ATMP Mycoplasma	25/100 tests
SMB95-1005/1006	Microsart® Research Mycoplasma	25/100 tests
SMB95-1007	Microsart® ATMP Sterile Release	10 samples
SMB95-1008	Microsart® ATMP Bacteria	100 tests
SMB95-1009	Microsart® Research Bacteria	25 tests
SMB95-1012	Microsart® ATMP Fungi	100 tests
SMB95-1014/1013	Microsart® Research Fungi	25/100 tests

Microsart® Calibration Reagent, 10⁸ genomes / vial, 1 vial (bacteria, including Mollicutes)

SMB95-2021	Mycoplasma arginini
SMB95-2022	Mycoplasma orale
SMB95-2023	Mycoplasma gallisepticum
SMB95-2024	Mycoplasma pneumoniae
SMB95-2025	Mycoplasma synoviae
SMB95-2026	Mycoplasma fermentans
SMB95-2027	Mycoplasma hyorhinis
SMB95-2028	Acholeplasma laidlawii
SMB95-2029	Spiroplasma citri
SMB95-2030	Bacillus subtilis
SMB95-2031	Pseudomonas aeruginosa
SMB95-2032	Kocuria rhizophila
SMB95-2033	Clostridium sporogenes
SMB95-2034	Bacteroides vulgatus
SMB95-2035	Staphylococcus aureus
SMB95-2036	Mycoplasma salivarium

Microsart® Calibration Reagent, 10⁶ genomes / vial, 1 vial (fungi)

SMB95-2044	Candida albicans
SMB95-2045	Aspergillus brasiliensis
SMB95-2046	Aspergillus fumigatus
SMB95-2047	Penicillium chrysogenum
SMB95-2048	Candida glabrata
SMB95-2049	Candida krusei
SMB95-2050	Candida tropicalis

Microsart® Validation Standard, 100 CFU / vial, 3 vials each (Mollicutes)

SMB95-2051	Mycoplasma orale
SMB95-2052	Mycoplasma pneumoniae

Microsart® Validation Standard, 99 CFU / vial, 6 vials each (bacteria* and fungi)

SMB95-2005	Bacillus subtilis
SMB95-2006	Pseudomonas aeruginosa
SMB95-2007	Kocuria rhizophila

SMB95-2008	Clostridium sporogenes
SMB95-2009	Bacteroides vulgatus
SMB95-2010	Staphylococcus aureus
SMB95-2037	Candida albicans
SMB95-2038	Aspergillus brasiliensis
SMB95-2039	Aspergillus fumigatus
SMB95-2040	Penicillium chrysogenum
SMB95-2041	Candida glabrata
SMB95-2042	Candida krusei
SMB95-2043	Candida tropicalis

* except for Mollicutes

DNA Extraction Kit

SMB95-2001	Microsart® ATMP Extraction (for bacteria and fungi)	50 extractions
SMB95-2003	Microsart® AMP Extraction (for mycoplasma)	50 extractions
SMB95-4000	Microsart® Proteinase K	50 extractions

Cleaning Spray

SMB95-5001	DNA Decontamination Reagent, spray bottle	250 ml
SMB95-5002	DNA Decontamination Reagent, refill canister	5 l

Cleaning Wipes

SMB95-5003	DNA Decontamination Reagent, wipes	50 wipes
SMB95-5004	DNA Decontamination Reagent, refill sachets	5 × 50 wipes

Limited Product Warranty

This warranty limits our liability for replacement of this product.

No warranties of any kind, express or implied, including, without limitation, implied warranties of merchantability or fitness for a particular purpose, are provided.

Trademarks

Microsart is a registered trademark of Sartorius.

Last technical revision: 2019-05-14

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Str. 20
37079 Goettingen, Germany

Phone +49 551 308 0
Fax +49 551 308 3289

 www.sartorius.com

Copyright by Sartorius, Goettingen, Germany.
All rights reserved. No part of this publication may be
reprinted or translated in any form or by any means
without the prior written permission of Sartorius.
The status of the information, specifications and
illustrations in this manual is indicated by the date
given below. Sartorius reserves the right to make
changes to the technology, features, specifications
and design of the equipment without notice.

Status:
September 2024,
Sartorius,
Goettingen, Germany

Printed in Germany on paper that
has been bleached without any use
of chlorine. | W
Publication No.: SM-6110-e240901
Order No.: 85037-546-75
Ver. 09 | 2024