

## Microsart® ATMP Extraction

Bacterial, fungi, and yeast DNA extraction kit  
Prod. No. SMB95-2001

Reagents for 50 reactions  
For use in research and quality control

Manufactured by:



Minerva Biolabs GmbH | Schkopauer Ring 13 | 12681 Berlin | Germany

## Symbols

---

**LOT**

Lot No.

---

**REF**

Order No.

---



Expiry date

---



Store at

---



Contains reagents for  
50 extractions

---



Manufacturer

---

# Contents

1. Intended Use.....	5
3. Test Principle.....	5
4. Notes on the Test Procedure.....	6
4.1 Handling and equipment recommendations.....	7
5. Reagents.....	8
6. Needed but not included.....	9
7. Specimen.....	12
Sample collection and storage.....	12
8. Test Procedure.....	13
8.1 Recommendation for product release testing.....	13
8.2 DNA extraction process.....	13
9. Related Products.....	14
Short Instructions.....	inside back cover



# 1. Intended Use

Microsart® ATMP Extraction kit introduces a state-of-the-art DNA extraction method for the isolation of bacterial, fungal, yeast DNA from cell cultures and cell culture-derived biologicals, like autologous transplants and other advanced therapy medicinal products (ATMP). The kit allows obtaining DNA templates for subsequent DNA amplification via PCR (e.g. for contamination detection, product release testing etc.).

# 2. Explanation of the Test

To achieve highest sensitivity and to avoid inhibitory effects in PCR testing, a DNA extraction step is recommended. For most test materials, DNA extraction methods, which provide suitable templates for PCR, are available. However, most of the DNA extraction kits available on the market are not free of DNA contaminations.

Microsart® ATMP Extraction introduces a unique DNA extraction method, which eliminates the risk of DNA contaminations, facilitating the detection of bacteria, fungi, and yeasts in cell culture and ATMPs via PCR.

The extraction procedure can be performed within 1 hour. In contrast to culture-based methods, for PCR tests, samples do not need to contain living material, as all intact particles (e.g. live, dormant, non-culturable etc.) are detected.

# 3. Test Principle

Microsart® ATMP Extraction kit was optimized for the extraction of genomic bacterial or fungal DNA from different sample matrices, including cell culture samples. The contamination risk has been minimized thanks to a reduced number of handling steps and a straightforward protocol.

An internal control DNA, provided with the Microsart® ATMP Bacteria or Microsart® ATMP Fungi kit, can be added to each sample prior to DNA extraction to monitor the extraction process. This will allow detection of false negative results, which may occur due to improper DNA extraction or PCR inhibition. Alternatively, the internal control DNA can be added directly to the master mix during PCR setup.

## 4. Notes on the Test Procedure

1. For *in vitro* use in research and quality control. This kit may be disposed of according to local regulations.
2. This kit should be used by trained staff, only. A clean lab coat and disposable gloves should be worn at all times while performing the assay.
3. To avoid DNA cross-contaminations, the complete test must be performed under sterile and DNA-free conditions (see chapter 4.1 for detailed information).
4. In case of work with living strains, the local regulation for S2 labs must be followed.
5. Aliquoting and repeatedly freezing and thawing the samples can increase the risk of sample contaminations. Therefore, this should be avoided whenever possible.
6. This extraction kit has been validated with 1 ml starting volume. If you use less than 1 ml it must be ensured that 99 CFU can be detected in the appropriate volume.
7. This kit is not validated for the extraction of mycoplasma DNA.
8. This leaflet must be widely understood for a successful use of Microsart® ATMP Extraction kit. The reagents supplied should not be mixed with reagents from different lots but used as an integral unit. The reagents of the kit should not be used beyond their shelf life.
9. Any deviation from the test method can affect the results.
10. For each test setup, at least one negative extraction control should be included. Positive controls facilitate the evaluation of the test.
11. The controls should be handled in the same manner as the samples.
12. Participation in external quality control programs, such as those offered by Minerva Biolabs GmbH ([www.minerva-biolabs.com](http://www.minerva-biolabs.com)), is recommended.

## 4.1 Handling and equipment recommendations

To avoid false positive results due to improper handling, the following actions are recommended:

1. To perform the test under sterile and DNA-free conditions, we recommend the use of an isolator/glovebox with an airlock.
2. The isolator/glovebox should be cleaned thoroughly with PCR Clean™ (Minerva Biolabs, Prod. No. 15-2025) or PCR Clean™ Wipes (Minerva Biolabs, Prod. No. 15-2001) before and during the working process.
3. All materials which are introduced into the isolator/glovebox should be cleaned thoroughly with PCR Clean™. Do not forget to clean the airlock with PCR Clean™. Pipettes and gloves should be cleaned thoroughly with PCR Clean™ Wipes prior to and during the process.
4. Avoid working above open tubes and avoid air turbulences due to rapid movements.
5. Be careful when opening the tubes. Do not touch the inner surface of the lid.
6. Always use a new, unopened DNA-free pipette filter tip-box for each assay. Reaction vials should be closed immediately after every pipetting step.

## 5. Reagents

Each kit contains reagents for 50 extractions. The expiry date of the unopened package is marked on the package label. The kit components are stored at ambient temperature until use. The lot specific Certificate of Analysis can be downloaded from the manufacturer's website ([www.minerva-biolabs.com](http://www.minerva-biolabs.com)).

<b>Kit Component Label Information</b>	<b>Quantity</b>	<b>Cap Color</b>
	<b>50 Extractions Order No. SMB95-2001</b>	
Lysis Buffer	2 × 13 ml	transparent
Suspension Buffer	4 × 1.5 ml	violet
Processing Tubes	50 × tubes	



## 6. Needed but not included

Microsart® ATMP Extraction kit contains reagents for sample collection and DNA extraction. General industrial supplies and reagents, usually available in PCR laboratories are not included:

### Consumables

- Laboratory gloves
- PCR Clean™ (Minerva Biolabs, Prod. No. 15-2025) and PCR Clean™ wipes (Minerva Biolabs, Prod. No. 15-2001)
- DNA-free pipette filter tips (we recommend Biosphere® filter tips from Sarstedt: 0.5-20 µl, Prod. No. 70.1116.210; 2-100 µl, Prod. No. 70.760.212; 20-300 µl, Prod. No. 70.765.210; 100-1000 µl, Prod. No. 70.762.211)
- DNA-free PCR reaction tubes (PCR 8-SoftStrips with attached caps from Biozym are recommended: 0.1 ml Low Profile, Prod. No. 710975 and 0.2 ml High Profile, Prod. No. 710970)

### Equipment

- Isolator/glovebox (further information, supplier and prices are available on request, please contact PCR@sartorius.com)
- Heat block with optional shaking function
- Microcentrifuge for 1.5 ml reaction tubes (Centrisart A-14, Prod. No. A-14-1EU)
- Vortex
- 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
- Rack for 1.5 ml tubes

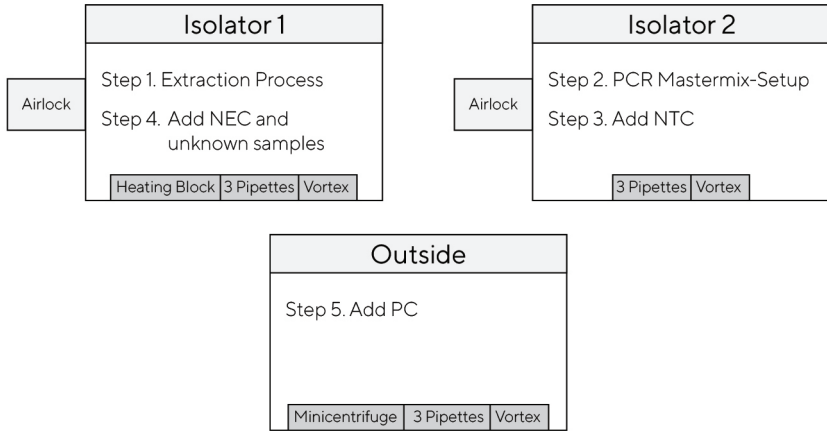
For subsequent PCR analysis, the following consumables and equipment are additionally required:

- PCR-based system for the detection of bacterial DNA: we recommend the Microsart® ATMP Bacteria kit (Sartorius Prod. No. SMB95-1008), or the Microsart® RESEARCH Bacteria kit (Sartorius Prod. No. SMB95-1009).
- PCR-based system for the detection of fungi and yeasts DNA: we recommend the Microsart® ATMP Fungi kit (Sartorius Prod. No. SMB95-1012), or the Microsart® RESEARCH Fungi kit (Sartorius Prod. No. SMB95-1013/1014).

**Note:** All these kits are qPCR-based assays, designed and optimized for the detection in cell cultures and cell culture-derived biologicals.

- qPCR device with filter sets for the detection of the fluorescent dyes FAM™ and ROX™ and suitable for 25 µl PCR reaction volumes.
- DNA-free PCR reaction tubes compatible with your specific qPCR device (PCR 8-SoftStrips with attached caps from Biozym are recommended: 0.1 ml Low Profile, Prod. No. 710975 and 0.2 ml High Profile, Prod. No. 710970).
- Minicentrifuge for PCR tubes
- Isolator/glovebox (for PCR setup)
- Vortex
- Set of 3 pipettes (see previous page)
- Rack for 1.5 ml tubes and for PCR tubes

Schematic overview of technical setup and experimental design:



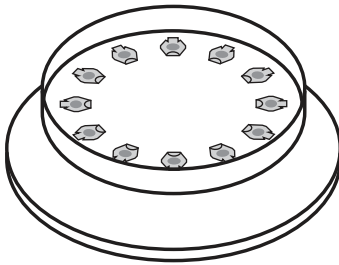
It is also possible to connect Isolator 1 and Isolator 2 via an airlock so that you can transfer the PCR tubes after Step 3 directly from Isolator 2 into Isolator 1. Please note that in this case you would need an additional airlock for Isolator 2.

# 7. Specimen

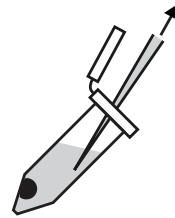
## Sample collection and storage

The kit has been validated using a maximum cell concentration of  $10^6$  cells/ml. Notably, the assay can be performed with different types of cell culture-derived material. Therefore, the optimal sampling parameters, like volume or cell number, can vary according to the specific characteristics of the sample (e.g. medium, cell type) and may require optimization of the procedure.

- 
1. max. 1 ml of cell culture or cell culture supernatant liquid material is transferred into a provided DNA-free 1.5 ml processing tube.
- 
2. Spin down for 15 min at a speed of at least  $16,200 \times g$  to sediment particles.  
**Attention:** Make sure to position the tubes in the centrifuge so to obtain a pellet on the back side of the tube, as described in the figure below (left).
- 
3. Carefully and completely discard the supernatant as described in the figure below (right). Proceed to DNA extraction. If DNA extraction cannot be performed immediately, freeze the samples at  $\leq -18 \text{ }^\circ\text{C}$ . Repeated freezing and thawing should be avoided.  
**Attention:** Samples can only be inactivated or frozen after this sample collection step.
- 



Make sure to position the tubes with the back side toward the outside of the rotor in order to obtain a pellet on the back side of the tube wall.



Slowly discard all the supernatant without disturbing the pellet.

## 8. Test Procedure

### 8.1 Recommendation for product release testing

The extraction process should be carried out with a negative extraction control (NEC) and samples in duplicates (= 3 extractions for 1 sample/product).

### 8.2 DNA extraction process

---

1. Add 500 µl Lysis Buffer (transparent cap) to cell pellet.

**Recommended for users of Microsart® ATMP Bacteria and Microsart® ATMP Fungi detection kit:**

The Internal Control DNA, which is included in the detection kit can also be used to monitor the extraction process.

Add 20 µl of Internal Control DNA to the sample, vortex briefly and proceed with step 2 as described. No additional Internal Control DNA is required for the PCR reaction mix.

---

2. Vortex vigorously for at least 30 sec until the pellet is completely dissolved.
- 

3. Heat at 80 °C (preferably with shaking) for 10 min.
- 

4. Spin down at 16,200 × g for 10 min.

**Attention:** Make sure to position the tubes in the rotor as indicated in the figure in chapter „Specimen“.

---

5. Remove supernatant carefully and completely following the instructions given in chapter „Specimen“. Make sure not to disturb or aspirate the pellet in the process.

**Attention:** There is a higher risk of PCR inhibition if liquid residues remain in the tube.

---

6. Add 100 µl Suspension Buffer (violet cap) and dissolve the DNA by thorough vortexing.
- 

Extracts can be stored for 6 days at +2 to +8 °C. If long-term storage is required, store at ≤ -18 °C. Repeated freezing and thawing should be avoided.

## 9. 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, 10<sup>8</sup> 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, 10<sup>6</sup> 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, 10 CFU / vial, 3 vials each (Mollicutes)

SMB95-2011	Mycoplasma arginini
SMB95-2012	Mycoplasma orale
SMB95-2013	Mycoplasma gallisepticum
SMB95-2014	Mycoplasma pneumoniae
SMB95-2015	Mycoplasma synoviae
SMB95-2016	Mycoplasma fermentans

SMB95-2017	Mycoplasma hyorhinis
SMB95-2018	Acholeplasma laidlawii
SMB95-2019	Spiroplasma citri
SMB95-2020	Mycoplasma salivarium

**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**

SMB95-2003	Microsart® AMP Extraction (for mycoplasma)	50 extractions
56-0002	Proteinase K**	50 extractions

**PCR Clean™ \*\***

15-2025	DNA Decontamination Reagent, spray bottle	250 ml
15-2200	DNA Decontamination Reagent, refill bottles	4 × 500 ml

**PCR Clean™ Wipes\*\***

15-2001	DNA Decontamination Reagent, Wipes	50 wipes
15-2002	DNA Decontamination Reagent, refill sachets	5 × 50 wipes

\*\* Distributed by Minerva Biolabs

Notes



## **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. Minerva Biolabs shall have no liability for any direct, indirect, consequential, or incidental damages arising out of the use, the results of use, or the inability to use this product.


## **Trademarks**

Microsart is a registered trademark of Sartorius Stedim Biotech GmbH. PCR Clean is a trademark of Minerva Biolabs GmbH.

Last technical revision: 2019-11-15

Sartorius Stedim Biotech GmbH  
August-Spindler-Str. 11  
37079 Goettingen, Germany

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

 [www.sartorius.com](http://www.sartorius.com)

Copyright by Sartorius Stedim Biotech GmbH,  
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 Stedim Biotech GmbH. The status of the information, specifications and illustrations in this manual is indicated by the date given below. Sartorius Stedim Biotech GmbH reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Status:  
October 2020,  
Sartorius Stedim Biotech  
GmbH, Goettingen, Germany

Printed in Germany on paper that  
has been bleached without any use  
of chlorine. | W  
Material No.: 1000054395  
Ver. 10 | 2020

## 1. Sample Collection



+ 1 ml  
sample material

Processing tube



15 min  
 $\geq 16,200 \times g$



Discard  
supernatant



Store at  $\leq -18 \text{ }^\circ\text{C}$

or proceed to  
DNA extraction

## 2. DNA Extraction



+ 500  $\mu\text{l}$   
Lysis Buffer  
(transparent cap)



$\geq 30$  sec vigorously



80  $^\circ\text{C}$ , 10 min



$\geq 16,200 \times g$ , 10 min



Remove supernatant  
carefully



+ 100  $\mu\text{l}$  Suspension  
Buffer (violet cap)



$\geq 30$  sec vigorously



DNA ready for PCR



incubate



vortex



centrifuge



add

storage at +18 - +25  $^\circ\text{C}$   
(ambient temperature)

Optional:

Add Internal Control from Microsart® ATMP Bacteria or Microsart® ATMP Fungi

Sartorius Stedim Biotech GmbH  
August-Spindler-Str.11  
37079 Goettingen, Germany

Phone +49 551 308 0

Fax +49 551 308 3289

 [www.sartorius.com](http://www.sartorius.com)

Copyright by Sartorius Stedim Biotech GmbH,  
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 Stedim Biotech GmbH. The status of the information, specifications and illustrations in this manual is indicated by the date given below. Sartorius Stedim Biotech GmbH reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Status:

October 2020,

Sartorius Stedim Biotech

GmbH, Goettingen, Germany

Printed in Germany on paper that  
has been bleached without any use

of chlorine. | W

Material No.: 1000054395

Ver. 10 | 2020