

Microsart® Calibration Reagent

For fungi and yeast species

Prod. No. SMB95-2044 *Candida albicans*
Prod. No. SMB95-2045 *Aspergillus brasiliensis*
Prod. No. SMB95-2046 *Aspergillus fumigatus*
Prod. No. SMB95-2047 *Penicillium chrysogenum*
Prod. No. SMB95-2048 *Candida glabrata*
Prod. No. SMB95-2049 *Candida krusei*
Prod. No. SMB95-2050 *Candida tropicalis*

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



Content

Contents

1. Intended Use.....	5
2. Explanation of the Product	5
3. Principle	6
4. Notes on the Procedure	6
5. Reagents	7
6. Needed but not included.....	8
7. Test Procedure	9
7.1 Rehydration of the reagents.....	9
7.2 Preparation of the dilutions.....	9
7.3 PCR	10
7.4 Evaluation.....	10
8. Related Products.....	11
Notes.....	13

1. Intended Use

Microsart® Calibration Reagent is titrated genomic DNA and can be used as amplification and sensitivity control for endpoint PCR (gel-based evaluation). If quantitative PCR is required, titrated genomic DNA can be serially diluted and used as PCR templates to create standard curves. The software will generate a standard curve from the known DNA dilutions and use it to determine the DNA concentration of the unknown samples.

2. Explanation of the Product

This product provides isolated genomic DNA of a specific species. An early passage strain of the specific microorganism is inoculated in a suitable culture medium and harvested at the end of the logarithmic growth phase by repeated washing and centrifugation. Following the extraction, the DNA concentration is quantified photometrically (OD_{260/280}) and with a double-stranded DNA (dsDNA)-specific ultrasensitive fluorescent dye, calibrated with weight reference standards and controlled by qPCR (compared to exactly quantified calibrator plasmids). The DNA concentration is adjusted with TE80 buffer.

The following species are available:

Candida albicans	ATCC 10231; DSM 1386
Candida glabrata	ATCC 90030; DSM 11226
Candida krusei	ATCC 6258; ATCC 749; DSM 6128; DSM 11956
Candida tropicalis	ATCC 4563; ATCC 7349; ATCC 450; DSM 11953
Aspergillus fumigatus	ATCC 9197; DSM 819
Aspergillus brasiliensis	ATCC 16404; DSM 1988
Penicillium chrysogenum	ATCC 9178; DSM 848

3. Principle

Each vial contains 1×10^6 copies of the complete fungal genome. The material cannot be used for cultivation methods. The genome standard can directly be used for PCR. Please note: For process validation alternatively the corresponding Microsart® Validation Standards can be used (see Related Products).

4. Notes on the Procedure

1. For in vitro use in research and quality control. This kit should be used only by trained persons. This kit does not contain hazardous substances and may be disposed of according to local regulations.
2. This leaflet must be widely understood for a successful use of the Microsart® Calibration Reagent. The supplied reagents should not be mixed with reagents from different lots and used as an integral unit. The reagents of the kit should not be used beyond their shelf life.
3. Any deviation from the described method can affect the results.
4. For each test setup, at least one negative control should be added.
5. Participation in external quality control programs, such as those offered by Minerva Biolabs GmbH (www.minerva-biolabs.com), is recommended.

5. Reagents

Each kit contains 1 vial of fungal DNA (1×10^6 genome copies) and 3 vials of buffer for the preparation of dilutions and negative controls. Samples are lyophilized for product stability reasons. The material has been inactivated prior to lyophilization and can be considered as non-infectious. The expiry date of the unopened product is specified on the package label. The kit components are stored at +2 to +8 °C upon arrival, at low humidity conditions. Lyophilized components must be stored ≤ -18 °C after rehydration.

Component Label Information	Order No.	Quantity	Cap Color
Calibration Reagent	SMB95-2044-	1 × lyophilized	green
Buffer	SMB95-2050	3 × 2.0 ml	white

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

6. Needed but not included

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

Consumables

- Laboratory gloves
- DNA-free pipette filter tips (Biosphere® filter tips from Sarstedt are recommended: 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)
- 1.5 ml reaction tubes, DNA- and RNA-free

Equipment

- Microcentrifuge for 1.5 ml reaction tubes
- 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

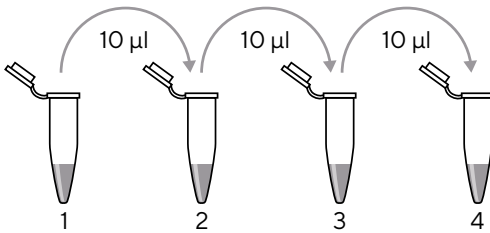
For PCR analysis, a fungi-specific DNA detection system is required. We recommend the Microsart® ATMP Fungi kit (Sartorius Prod. No. SMB95-1012), or the Microsart® Research Fungi kit (Sartorius Prod. No. SMB95-1013/1014).

7. Test Procedure

7.1 Rehydration of the reagents

All reagents must be equilibrated to room temperature prior to use.

1. Centrifuge the tube briefly to collect the lyophilized material at the bottom of the tube.
2. Add 100 μl Buffer (white cap) to the vial containing the Calibration Reagent (green cap) to obtain a concentration of 1×10^4 genomes/ μl .
3. Incubate 5 min at room temperature.
4. Vortex for 10 sec and spin for 5 sec with the “pulse” option or at $5000 \times g$.
5. Aliquot the vial content in DNA-free tubes and freeze $\leq -18^\circ\text{C}$ or directly proceed to step 7.2.



7.2 Preparation of the dilutions

1. Thaw the resuspended Calibration Reagent if necessary.
2. Label five 1.5 ml reaction tubes consecutively and fill each with 90 μl of the provided Buffer (white cap).
3. Vortex Calibration Reagent for at least 5 sec at medium speed.
4. Add 10 μl of the Calibration Reagent to reaction tube no. 1, close the tube and vortex briefly at medium speed.
5. Add 10 μl of the content of reaction tube no. 1 to reaction tube no. 2.
6. Close the tube and vortex briefly at medium speed.
7. Continue with the following reaction tubes of the dilution series as described above.

7.3 PCR

Please follow the instructions given in the manual of the specific PCR kit. The volume used as template for PCR defines the number of genome copies per reaction:

Reaction tube	2 µl sample volume	10 µl sample volume	50 µl sample volume
1.	2×10^3 genome copies	1×10^4 genome copies	5×10^4 genome copies
2.	200 genome copies	1000 genome copies	5000 genome copies
3.	20 genome copies	100 genome copies	500 genome copies
4.	2 genome copies	10 genome copies	50 genome copies
5.	-	1 genome copies	5 genome copies

7.4 Evaluation

In qPCR, the Ct values should decrease linearly with increasing DNA concentrations per reaction, when a suitable PCR assay is used. The software of the qPCR device builds a standard curve and calculates the slope using the DNA concentrations provided by the user and the obtained Ct values. Also, the unknown sample DNA concentrations are automatically determined by interpolation of the Ct values, obtained for these samples, on the standard curve. The following results represent an example of such evaluation and were generated using the CFX96 Touch™ instrument.

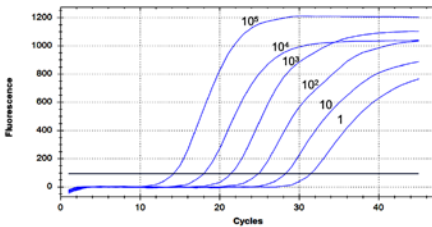


Fig. 1: Amplification curves of a dilution series from 10^5 to 1 genome copies/ reaction.

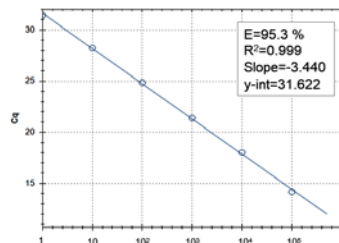


Fig. 2: Standard curve generated with CFX96 Touch™ instrument using second derivative maximum method and the data from Fig. 1.

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® 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

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

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

CFX96 Touch is a trademark of Bio-Rad Laboratories, Inc. Microsart is a registered trademark of Sartorius.

Last technical revision: 2019-11-19

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
DIR No.: 2543685-000-00
Publication No.: SM-2044-e240901
Ver. 09 | 2024