



sartorius

# Mycoplasma Assays for qPCR

## The new Gold Standard

Mycoplasma qPCR assays offer gold standard performance, design and configuration flexibility, affordability, and manufacturing quality.

### Flexibility

Select the product you need and take advantage of the protocol which fits perfectly to your sample simply by design.

- ▶ Fast and reliable screening of cell cultures in research, ATMPs or high end confidence testing in production – well designed kits for any need.
- ▶ Protocols for testing 2 µl, 200 µl or up to 18 ml are available.
- ▶ Open system, no hardware bundle – take advantage of the qPCR you already own and keep your investment low.

### Convenience

All three kits are based on easy-to-use real-time PCR technology for fast and reliable results.

- ▶ The TaqMan® assay design reduces the number of reactions needed, as it enables the use of internal controls. This increases the reliability and reduces the costs of the tests.
- ▶ The lyophilized Mycoplasma Mix contains all reaction components including the polymerase and can be easily reconstituted with the rehydration buffer included in the kit. It comes in aliquots of 25 reactions for highest convenience and long term stability.

### Performance

By teaming up with Minerva Biolabs, Sartorius has gained a partner with long standing experience and an extensive research background in qPCR – providing performance-guaranteed reagents and results.

- ▶ Comprehensive validation according to European Pharmacopoeia 2.6.7.
- ▶ Challenged in proficiency tests and in-depth robustness studies.



# Kit Selection

Features	Microsart® RESEARCH Mycoplasma	Microsart® ATMP Mycoplasma	Microsart® AMP Mycoplasma
Recommended use   scope	Testing of cell culture materials in research and development	Testing of autologous cell transplants (ATMPs)	Regulated in-process and lot-release testing
Type of PCR	5' nuclease assay for qPCR	5' nuclease assay for qPCR	5' nuclease assay for qPCR
Device requirements	None. Kit can be applied on any qPCR cycler suitable to detect FAM™ and ROX™ dyes	None. Kit can be applied on any qPCR cycler suitable to detect FAM™ and ROX™ dyes	Kit can be applied on any qPCR cycler suitable to detect FAM™ and ROX™ dyes and accepting 100 µl reaction volume
Kit components	<ul style="list-style-type: none"> <li>- lyophilized primers   nucleotides   probes   polymerase   internal amplification control, aliquoted in 25 reactions</li> <li>- rehydration buffer</li> <li>- lyophilized positive control</li> <li>- PCR grade water</li> </ul>	<ul style="list-style-type: none"> <li>- lyophilized primers   nucleotides   probes   polymerase, aliquoted in 25 reactions</li> <li>- internal amplification control usable as DNA extraction control</li> <li>- rehydration buffer</li> <li>- lyophilized positive control</li> <li>- PCR grade water</li> </ul>	<ul style="list-style-type: none"> <li>- lyophilized primers   nucleotides   probes   polymerase, aliquoted in 25 reactions</li> <li>- internal amplification control usable as DNA extraction control, aliquoted in 25 reactions</li> <li>- rehydration buffer</li> <li>- lyophilized positive control, aliquoted in 25 reactions</li> <li>- PCR grade water</li> </ul>
Package sizes	Cat.-No. SMB95-1005 25 tests Cat.-No. SMB95-1006 100 tests	Cat.-No. SMB95-1003 25 tests Cat.-No. SMB95-1004 100 tests	Cat.-No. SMB95-1001 25 tests Cat.-No. SMB95-1002 100 tests
Sample volume	2 µl	200 µl	200 µl to 18 ml
Sample volume per PCR	2 µl	10 µl	50 µl
EP 2.6.7 compliance	No	Yes, for ATMPs only	Yes, comprehensive
Validation	Specificity and sensitivity based on genome units	Specificity and sensitivity based on CFU for all EP listed mycoplasma species, robustness for ATMPs only	Comprehensive specificity testing including production cell lines and various bacteria species; sensitivity based on CFU for all EP listed mycoplasma species; robustness for different types of cell culture materials and supplements
Result evaluation	Cycler based, real-time PCR; validated protocols available	Cycler based, real-time PCR; validated protocols available	Cycler based, real-time PCR; validated protocols available
Required consumables	PCR reaction tubes	PCR reaction tubes	PCR reaction tubes
Optional consumables	None	<ul style="list-style-type: none"> <li>- 10CFU™ Sensitivity Standards for validation and performance control (e.g. SMB95-2011)</li> <li>- Microsart® AMP Extraction kit (Cat. No. SMB95-2003)</li> </ul>	<ul style="list-style-type: none"> <li>- 10CFU™ Sensitivity Standards for validation and performance control (e.g. SMB95-2011)</li> <li>- Microsart® AMP Extraction kit (Cat. No. SMB95-2003)</li> <li>- Vivaspin® 6 or 20 (e.g. VS0641, VS2041) in combination with Microsart® AMP Coating Buffer (SMB95-2002) to process volumes higher than 200 µl</li> </ul>
Shelf life and storage	Components can be stored at +2 to +8 °C for at least 12 months. After rehydration the reagents must be stored at -18 °C	Components can be stored at +2 to +8 °C for at least 12 months. After rehydration the reagents must be stored at -18 °C	Components can be stored at +2 to +8 °C for at least 12 months. After rehydration the reagents must be stored at -18 °C

FAM and ROX are trademarks of Applied Biosystems Corporation or its subsidiaries in the US and | or certain other countries.

Microsart® RESEARCH Mycoplasma, Microsart® ATMP Mycoplasma and Microsart® AMP Mycoplasma are distributed by Sartorius Lab Products and Services ([www.sartorius.com](http://www.sartorius.com)). Please contact your local Sartorius representative for local availability and ordering. Contact Sartorius Stedim Biotech or Minerva Biolabs for technical requests or any assistance.

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

Phone +49.551.308.0  
Fax +49.551.308.3289

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