

# SARTORIUS

## Simplifying Progress

### Microsart® Mycoplasma Kits

2022



# Microsart® Mycoplasma Kits



Features	Microsart® Research Mycoplasma	Microsart® ATMP Mycoplasma	Microsart® AMP Mycoplasma
Recommended use	Testing of cell culture materials in research and development	Testing of autologous cell transplants (ATMPs)	Regulated in-process and lot-release testing of cell cultures and cell-culture derived biologicals, with increased sample volume input
Type of PCR	5' nuclease assay for qPCR	5' nuclease assay for qPCR	5' nuclease assay for qPCR
Device requirements	Kit can be applied on any qPCR cyclers suitable to detect FAM™ and ROX™ dyes	Kit can be applied on any qPCR cyclers suitable to detect FAM™ and ROX™ dyes	Kit can be applied on any qPCR cyclers 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</li> <li>Rehydration buffer</li> <li>Lyophilized positive control</li> <li>PCR grade water</li> </ul>
Package sizes	Cat. No. SMB95-1005 25 tests	Cat. No. SMB95-1003 25 tests	Cat. No. SMB95-1001 25 tests
Sample volume	2 µL (optional 200 µL when combined with prior DNA extraction)	200 µL	200 µL to 18 mL
Sample volume per PCR	2 µL	10 µL	50 µL

**SARTORIUS**

**Mycoplasma Assays for PCR**  
**The New Gold Standard**

Mycoplasma PCR Assays Offer Gold Standard Performance, Design & Configuration Flexibility, and Quality

**Benefits**

- Select the product you need and take advantage of the protocol that 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 PCR instrument you already own and keep your investment low.
- Compatibility with many commercial qPCR machines with validated protocols.

# SARTORIUS

## Simplifying Progress

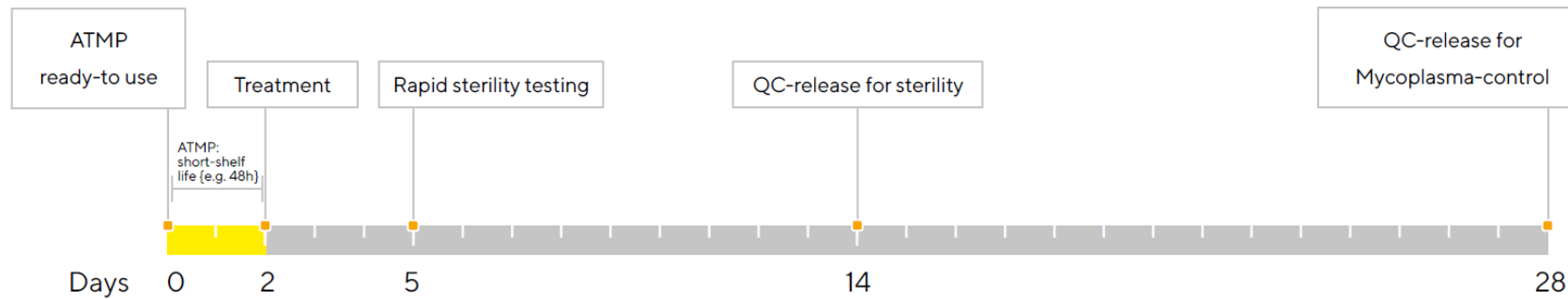
Microsart® AMP Extraction  
Microsart® ATMP Mycoplasma

2022

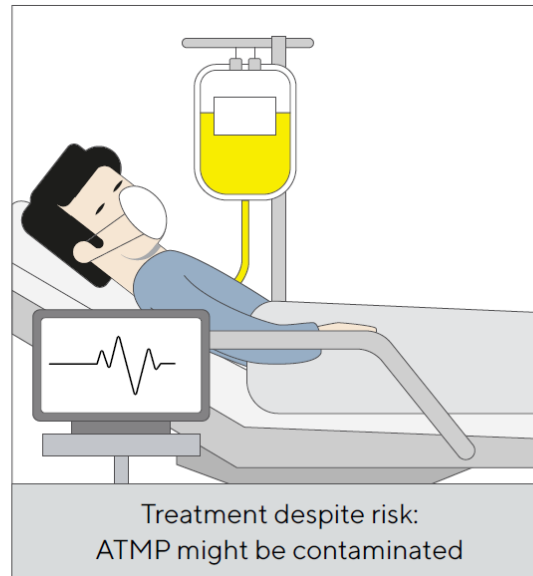
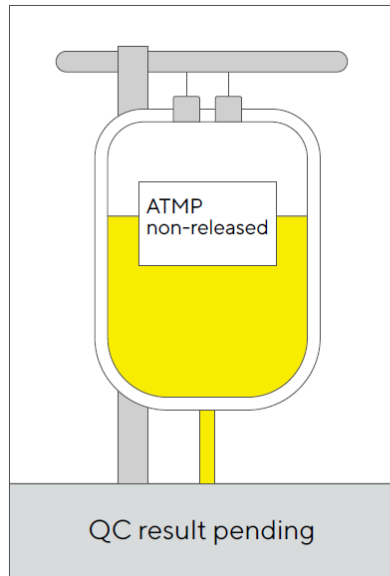


Why new methods?

# ATMPs put microbiological QC to novel challenges

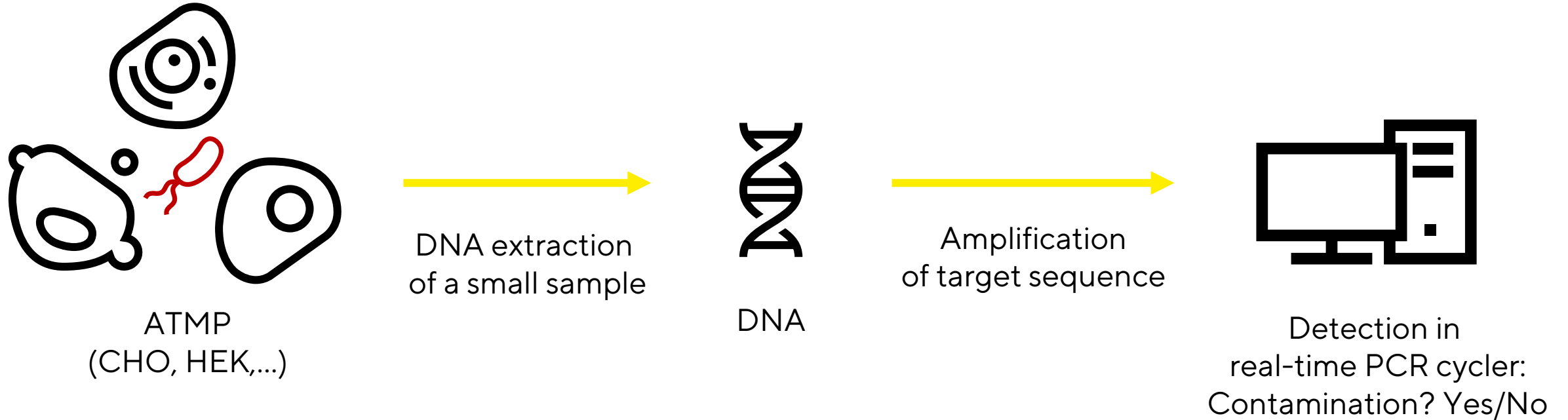


**5, 14 or 28 days  
of waiting  
is too long  
for ATMPs!**



Why new methods?

## Nucleic acid techniques



Results within 3 hours!

# Mycoplasma contamination detection

- Real-time PCR allows detection of Mycoplasma
  - In 3 h
  - Down to 5-10 CFU/ml
- Validated combination
  - In accordance with EP 2.6.7 and USP 63
- Support
  - **Product Validation Report** containing all experimental details
  - **Matrix Validation Proposal** giving an overview of the required set up and materials
  - **Matrix Validation Template** containing detailed information for the customer specific matrix validation
  - **Technical support** during matrix validation process



Microsart® AMP Extraction



Microsart® ATMP Mycoplasma

# Workflow Mycoplasma contamination detection

- DNA isolation using the column-based Microsart® AMP Extraction kit
  - Columns allow to isolate DNA from the whole ATMP sample (supernatant & cells)
  
- Real-time PCR using the Microsart® ATMP Mycoplasma kit
  - Taq-Man® System → reduce false-positive signals
  - Duplex assay → reduce false-negative signals
  - Universal assay for different real-time PCR cyclers → FAM™ and ROX™
  - High stability & no freezing → Lyophilized reagents



Microsart® AMP Extraction



Microsart® ATMP Mycoplasma



Simplifying Progress

Technical background  
DNA-based detection methods

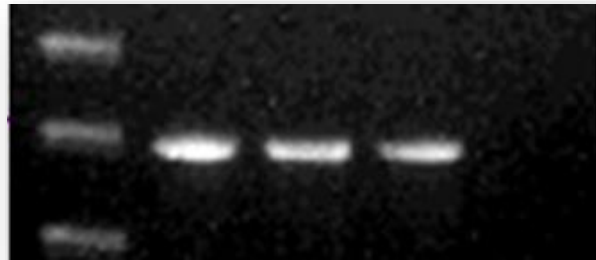
**SARTORIUS**



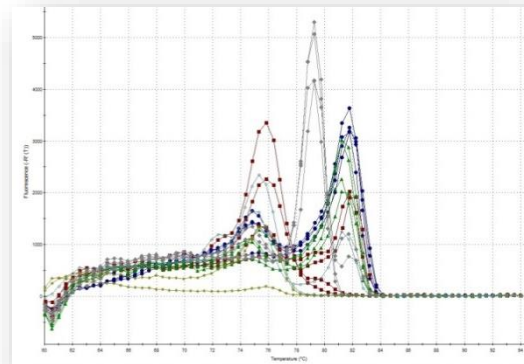
# DNA-based detection methods

real-time PCR

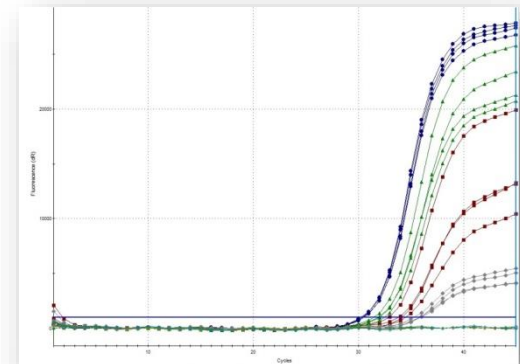
Conventional PCR



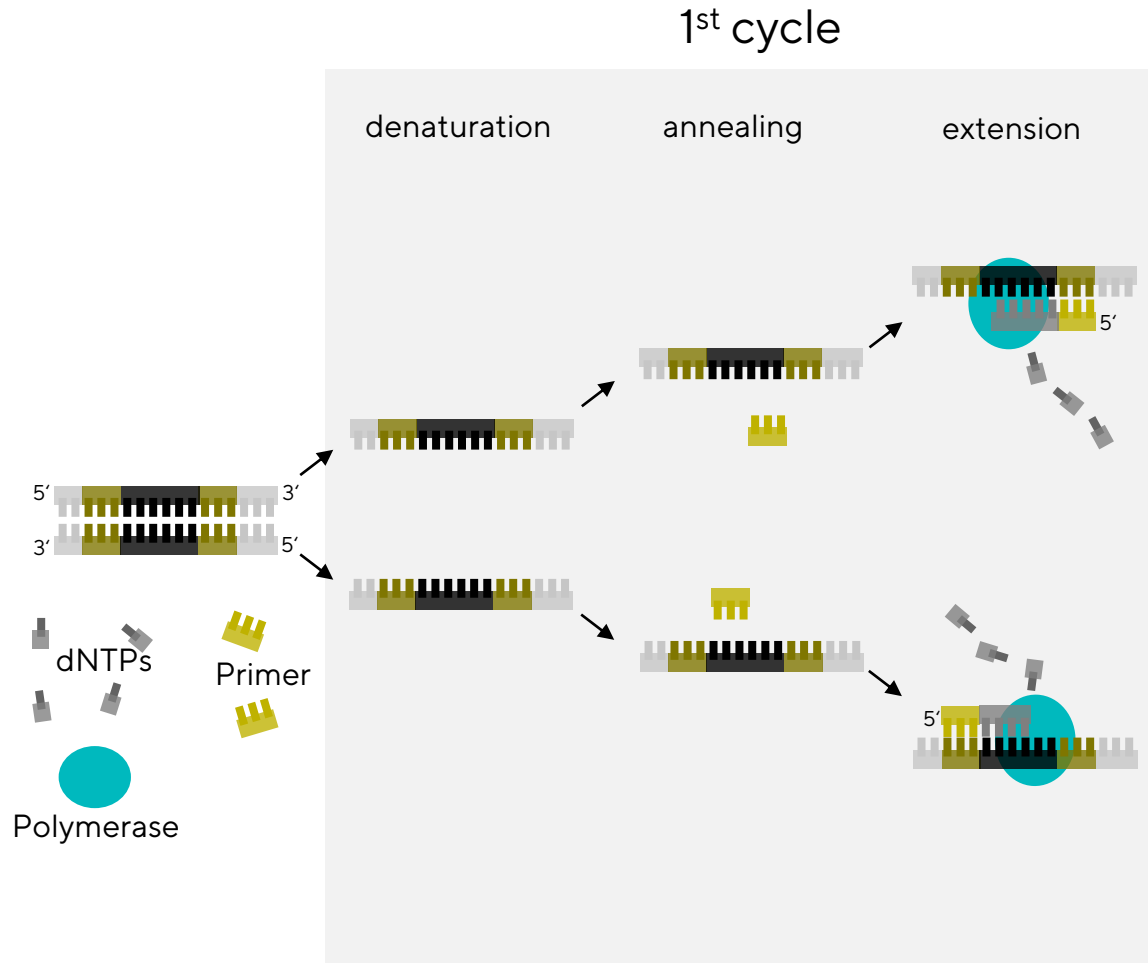
SYBR Green I



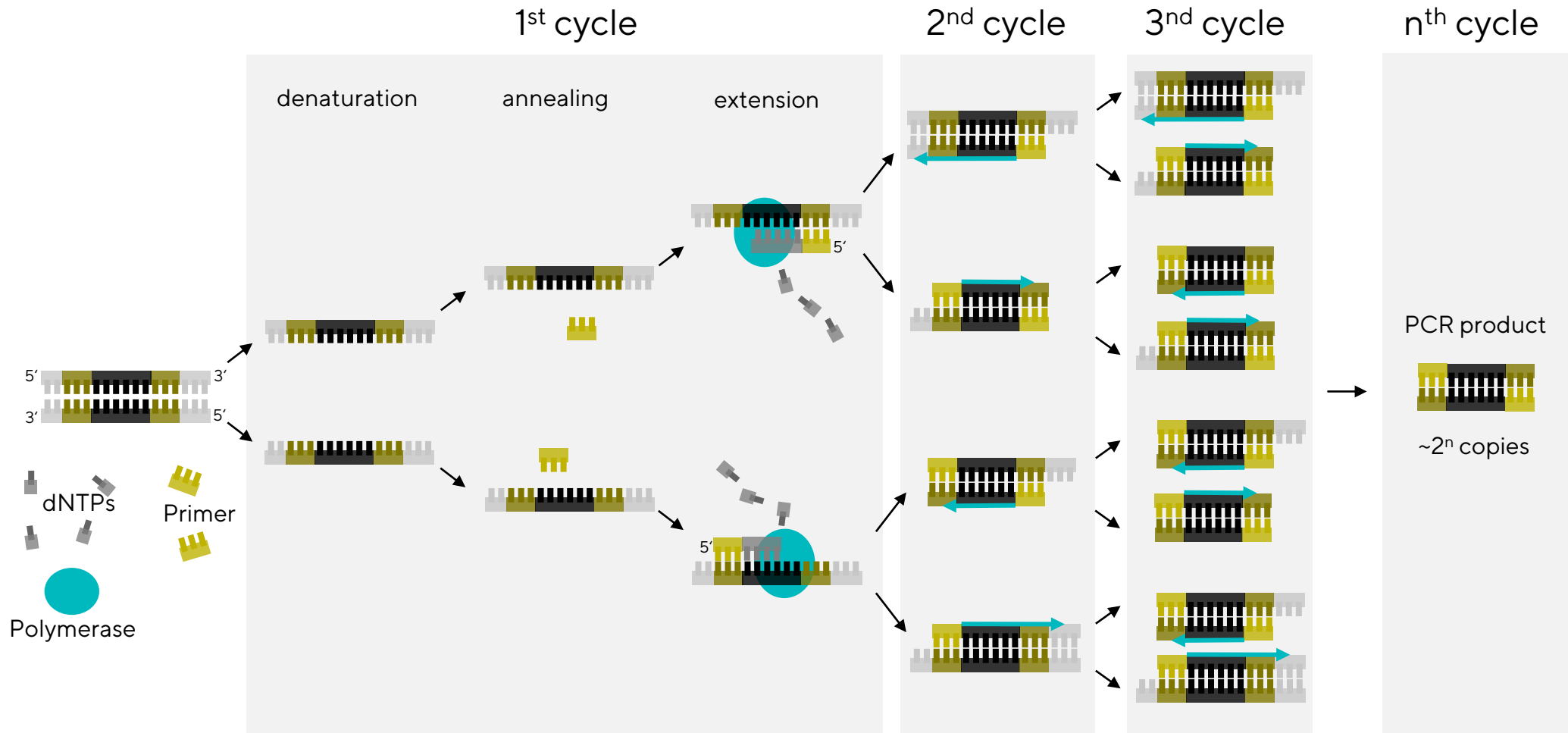
TaqMan® Probe



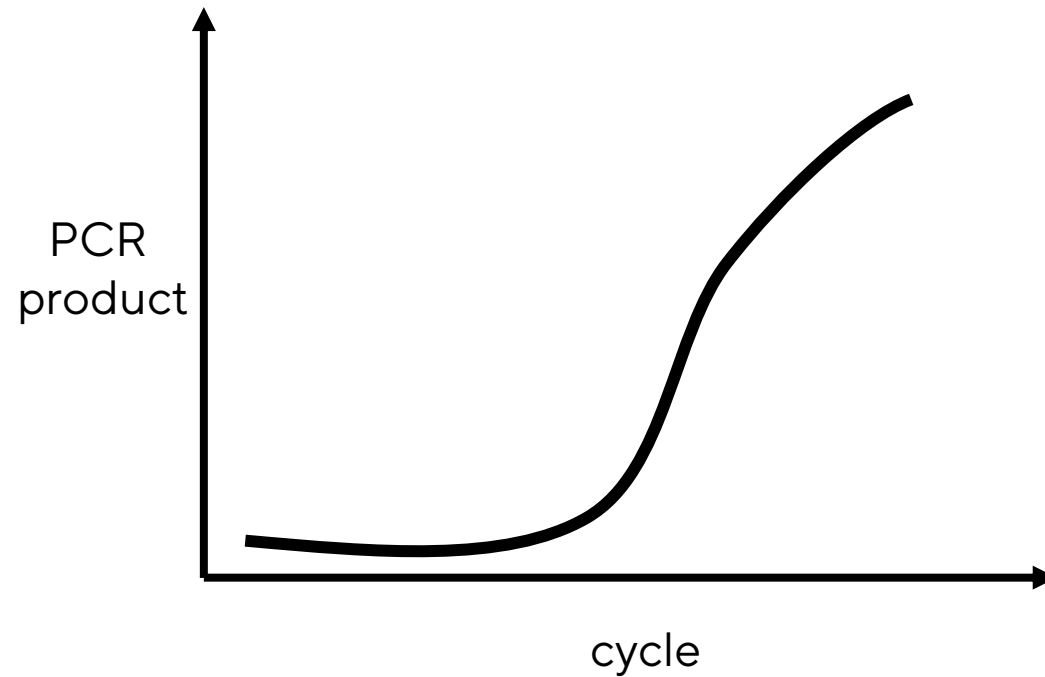
# What is a conventional PCR?



# What is a conventional PCR?

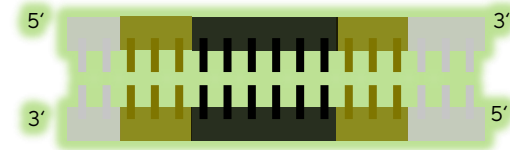
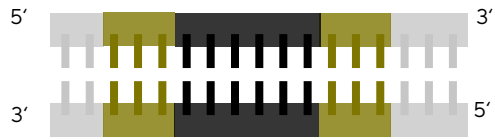
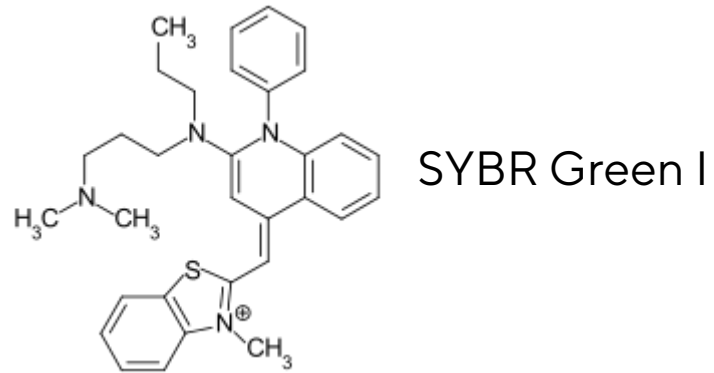


# A real-time PCR visualizes the reaction

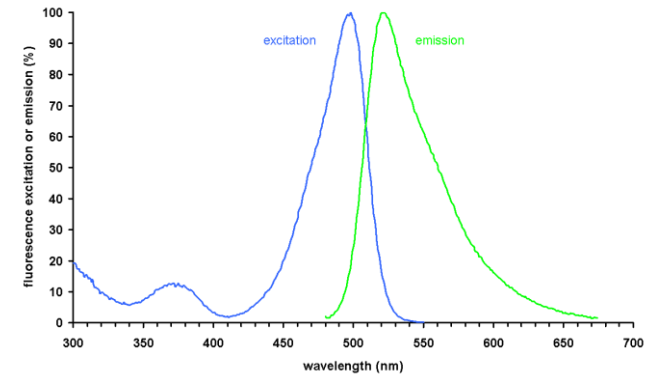


How does that work?

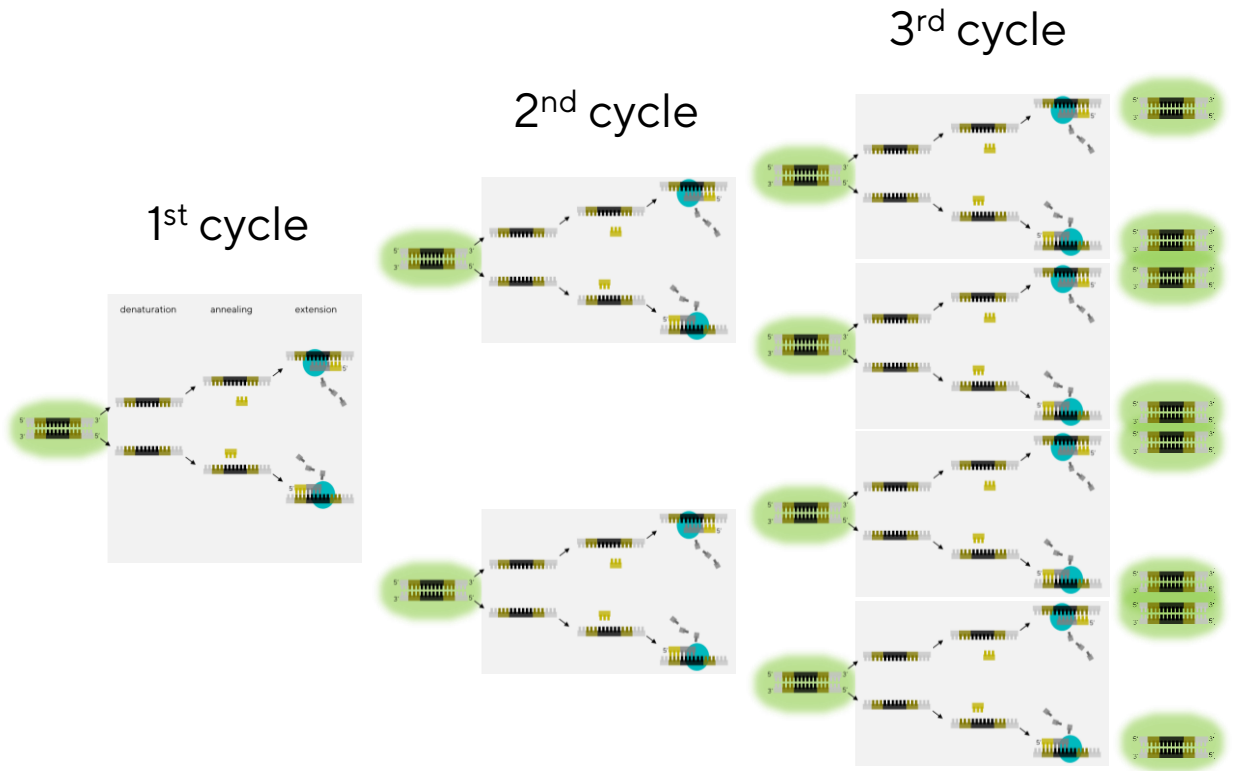
# Real-time PCR



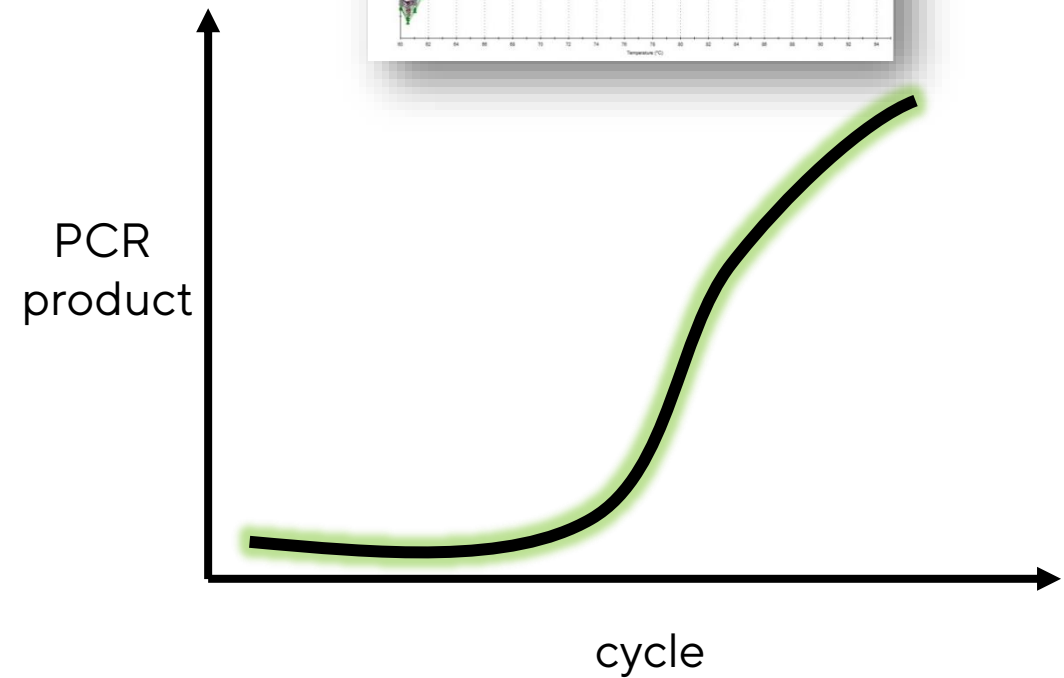
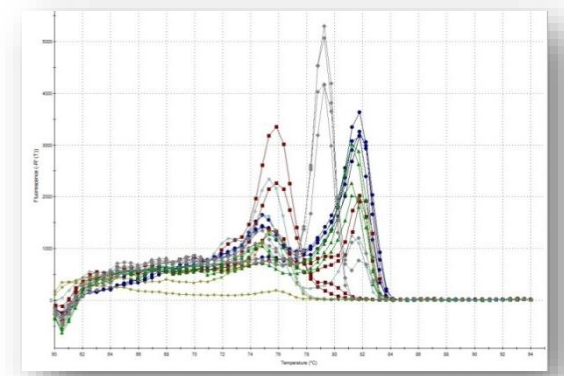
The dye SYBR Green I binds to double stranded DNA!



# Real-time PCR

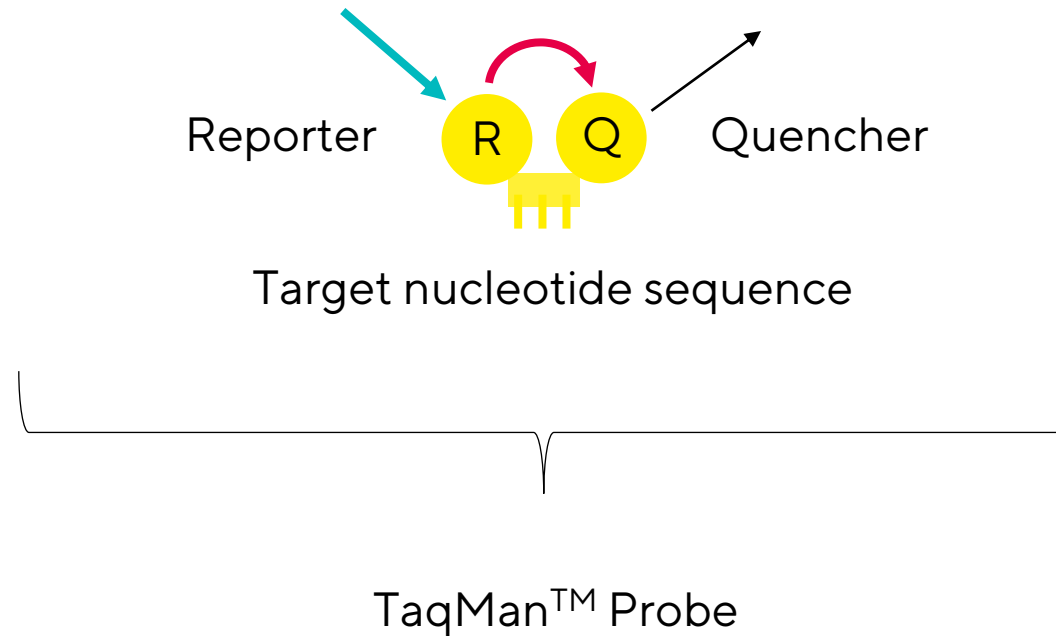


SYBR Green I



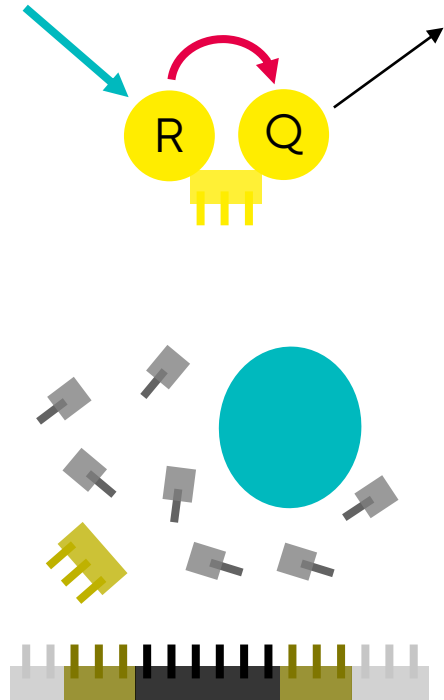
Unspecific binding of SYBR Green I can result in false-positive signals!

# A TaqMan™ probe is more specific compared to SYBR Green I

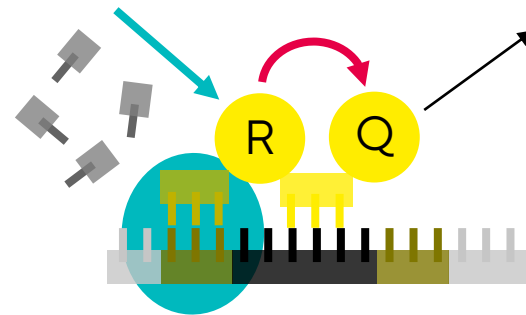


# TaqMan™ real-time PCR

TaqMan™ probe is degraded during real-time PCR

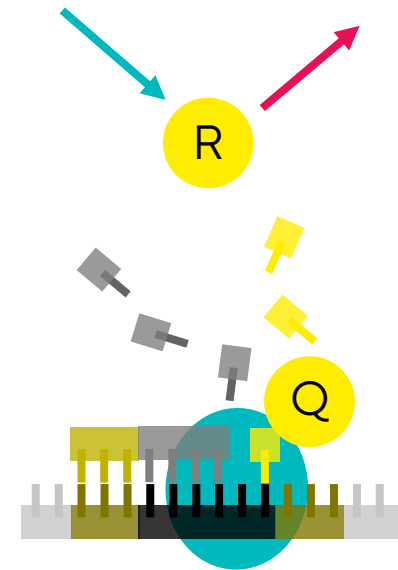


As long as the probe is complete  
no light signal can be detected



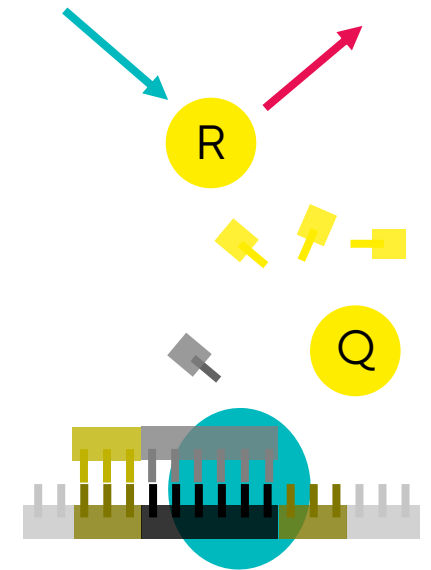
Taq Polymerase functions:

- DNA amplification
- 5'-3' exonuclease activity



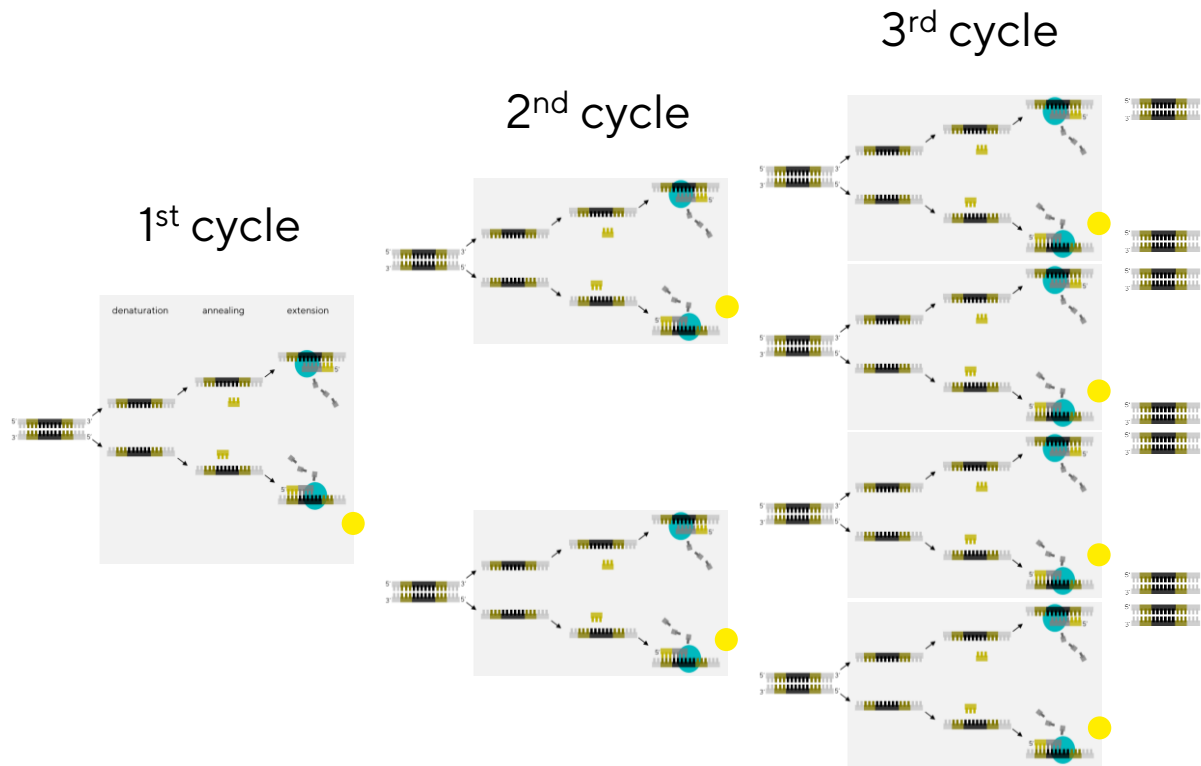
During elongation:

- Polymerase hydrolyses probe
- Dye and quencher are separated
- Reporter dye emits light signal

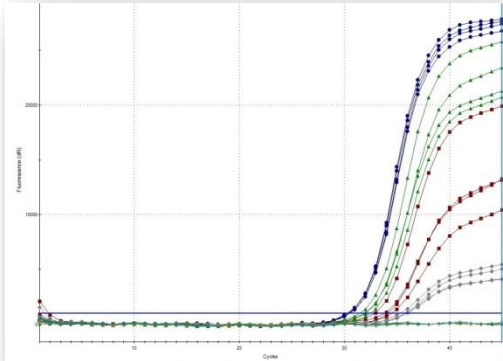




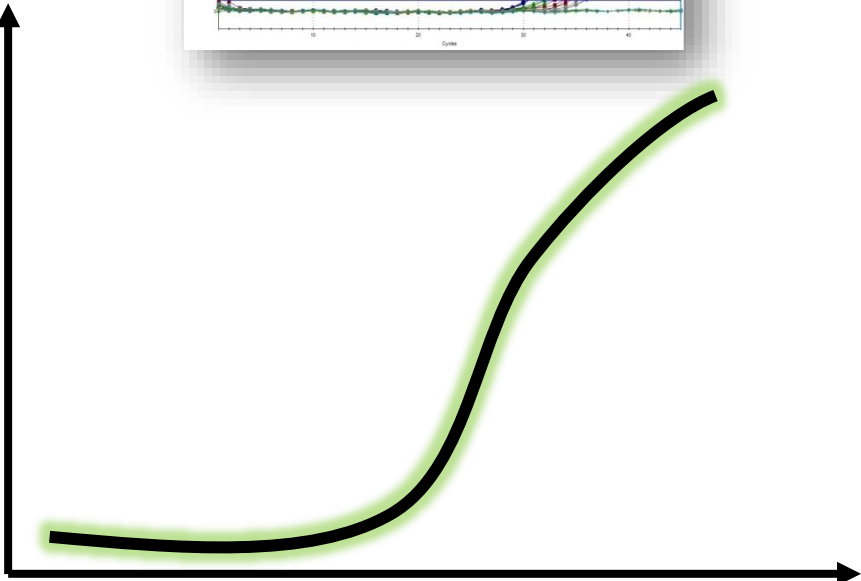
# TaqMan™ real-time PCR



TaqMan® Probe



PCR product



cycle

The specificity of TaqMan™ system **reduces false-positive results!**

# A duplex real-time PCR assay monitors PCR functionality

Problem:

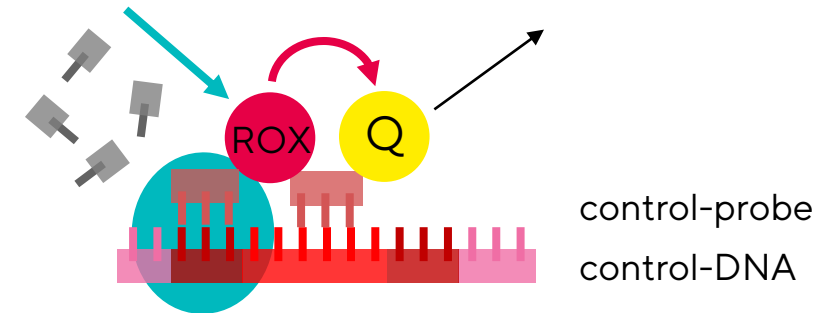
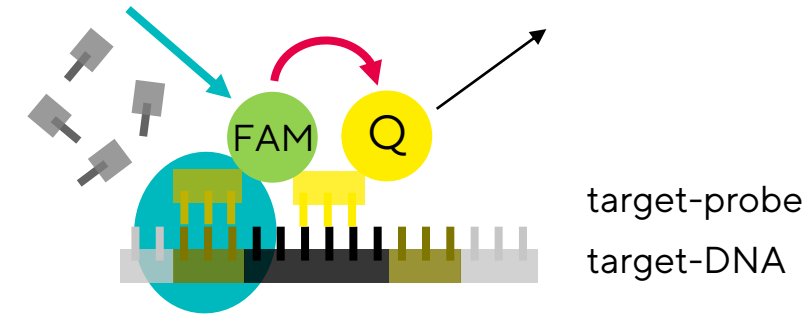
What does **no signal** mean?  
No microbial contamination?  
PCR inhibition?

Solution:

Include a second real-time PCR and  
a control DNA that must lead to a  
signal!

→ If this internal control reaction  
does not lead to a signal,  
the PCR is inhibited.

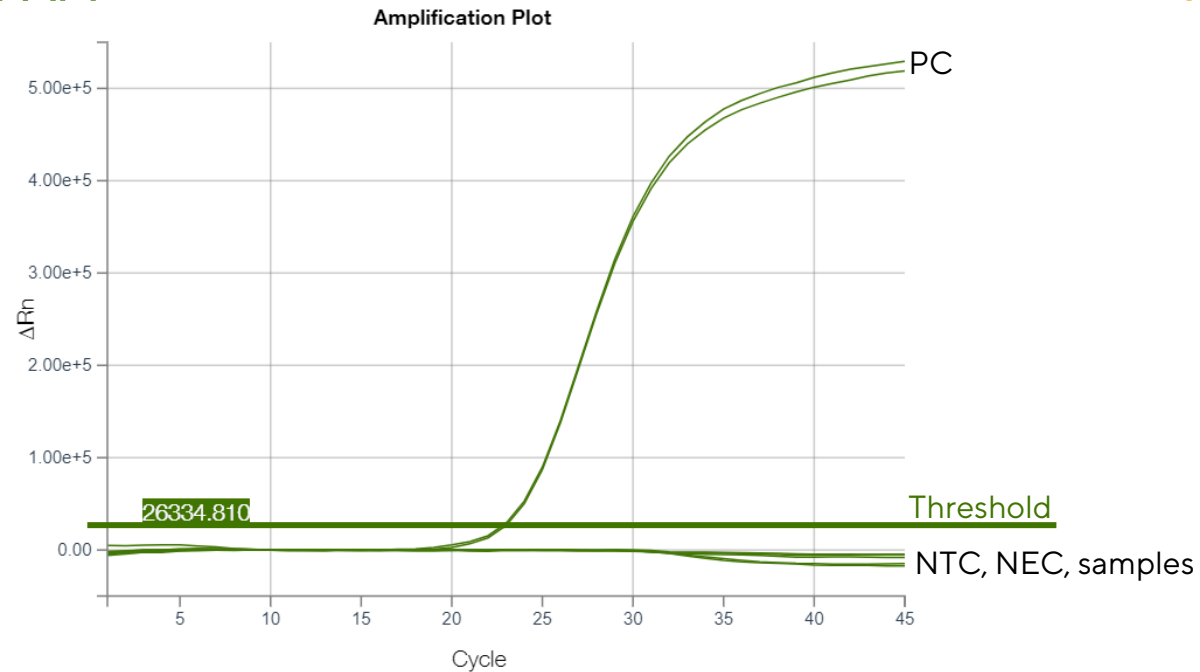
**Duplex assay**  
= two independent real-time PCRs  
in one run using different  
fluorophores



The internal control reaction **reduces false-negative results!**

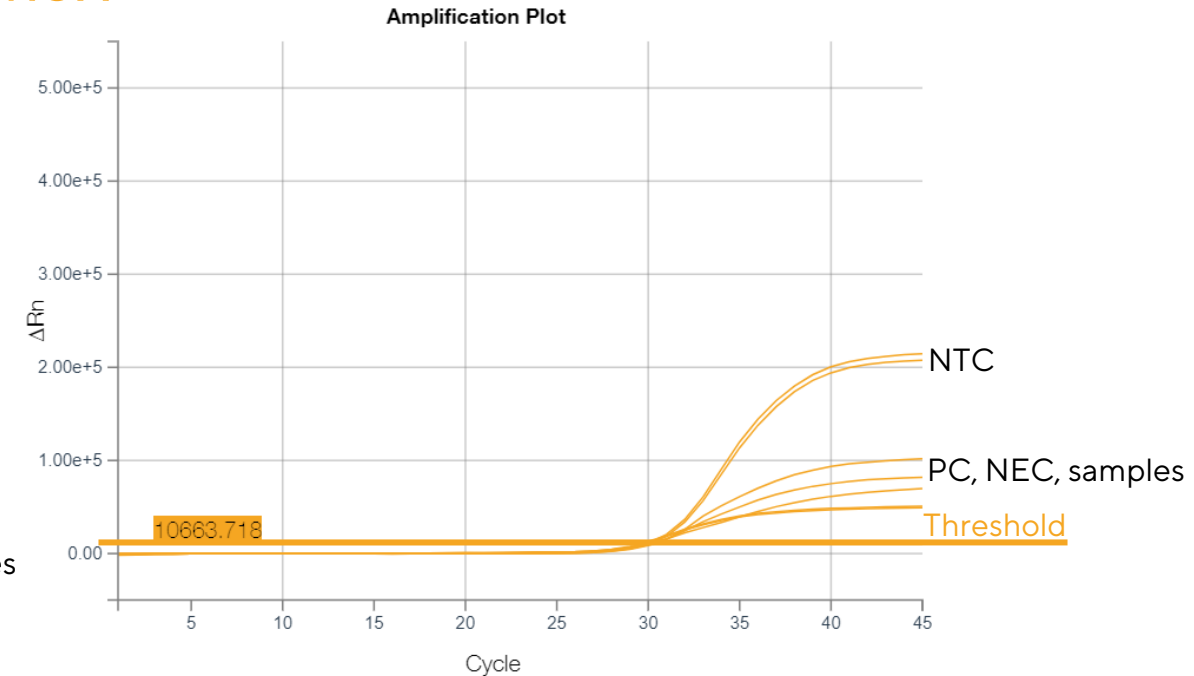
# A duplex real-time PCR Analysis

FAM



There is no contamination in the samples, because only the positive control is detectable in the FAM channel

ROX



There is no PCR inhibition, because the internal control DNA was detected in all reactions.

## What is in the kits?

### real-time PCR master mix

- Primer for target DNA
- Primer for control DNA
- FAM probe for target DNA
- ROX probe for control DNA
- Taq polymerase
- Buffer

Rehydration buffer

Internal Control DNA

Positive Control DNA

Ultrapure Water



Simplifying Progress

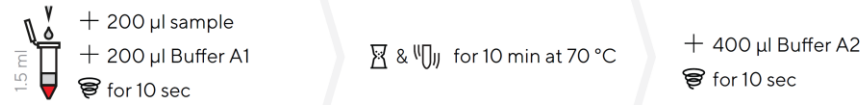


Hands on Mycoplasma detection

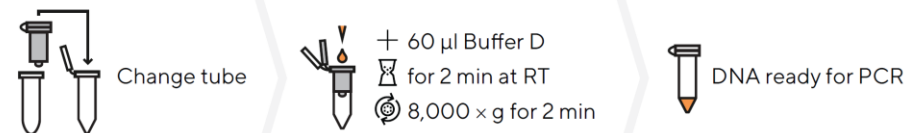
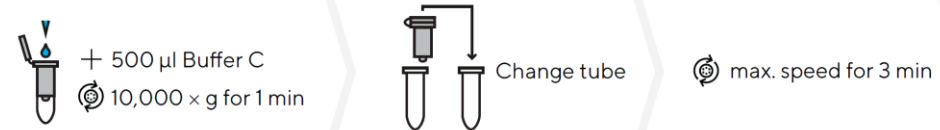
**SARTORIUS**

# Microsart® AMP Extraction

## 1. Preparation of Cell Culture Material

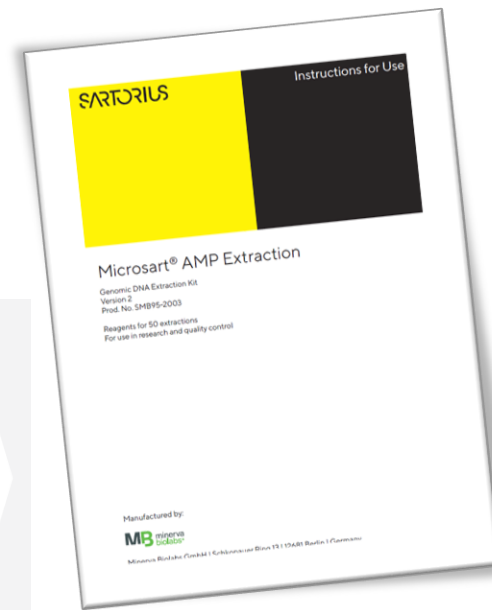


## 2. DNA Isolation



- ⌚ incubate
  - 🌀 vortex
  - 🌀 centrifuge
  - + add
  - 🌀 shake
- storage +18 - +25 °C (RT)

This procedure overview is not a substitute for the detailed manual.

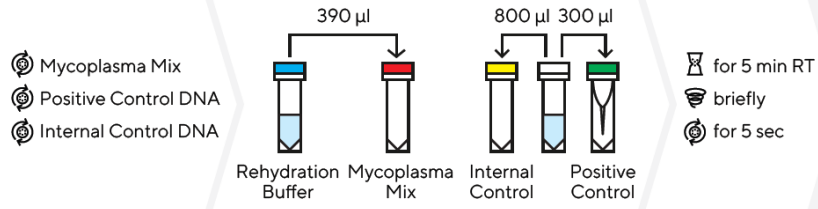


Silica membrane based extraction

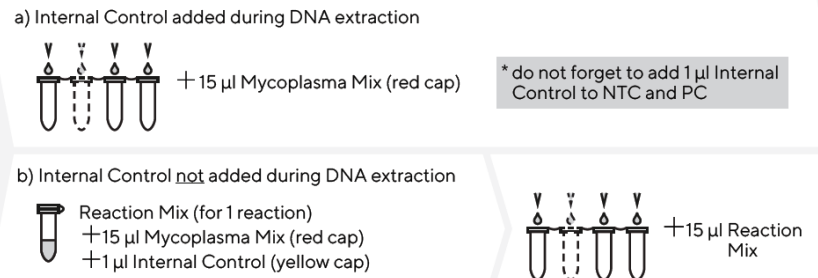
# Microsart® ATMP Mycoplasma



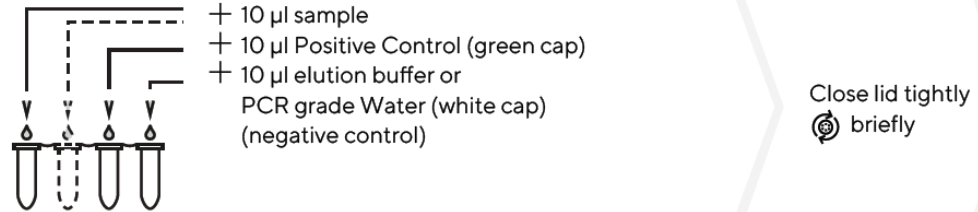
## 1. Rehydration of Reagents



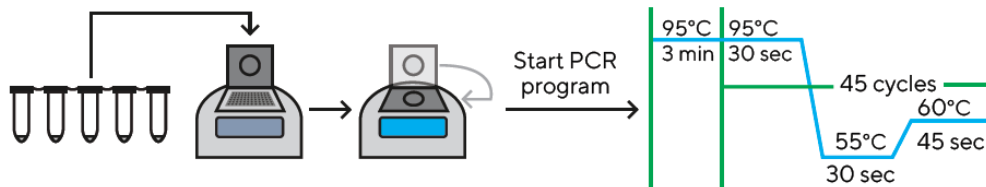
## 2. Preparation of PCR Reaction Mix



## 3. Addition of Samples and Controls



## 4. Start of the qPCR Reaction

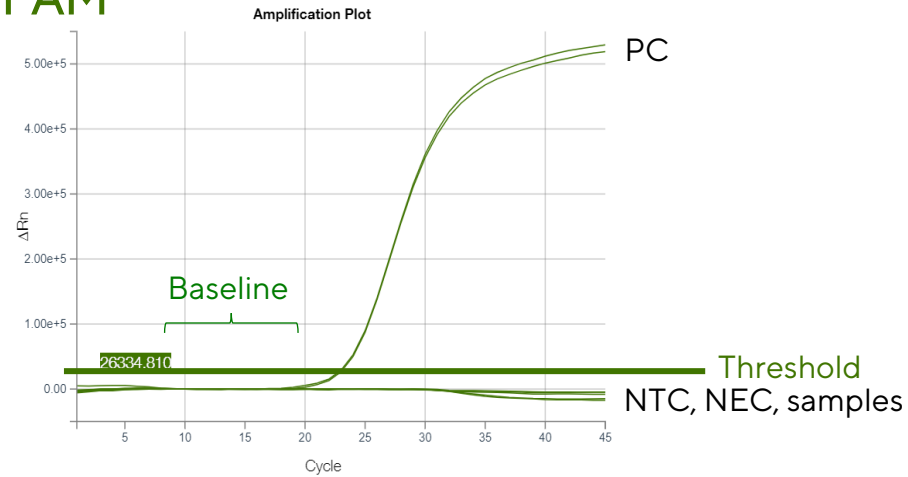


- Rehydration Buffer
- Mycoplasma Mix
- PCR grade Water
- Positive Control
- Internal Control
- ⏰ incubate
- 🌀 vortex
- 🌀 centrifuge
- ⊕ add
- storage +2 - +8 °C
- after rehydration ≤ -18 °C

This procedure overview is not a substitute for the detailed manual. ST\_SI\_Microsart®-ATMP-Mycoplasma\_03\_EN

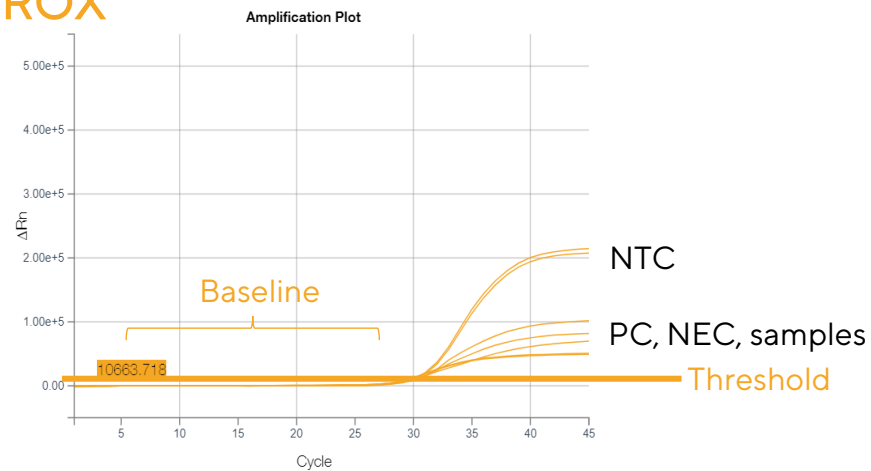
# Microsart® ATMP Mycoplasma - Analysis

FAM





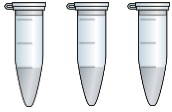


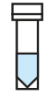












1. Set the baseline to level the curves
2. Set the threshold
3. Check if all controls are as expected  
→ see next slide
4. Analyze your samples

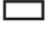


ROX





# Result interpretation

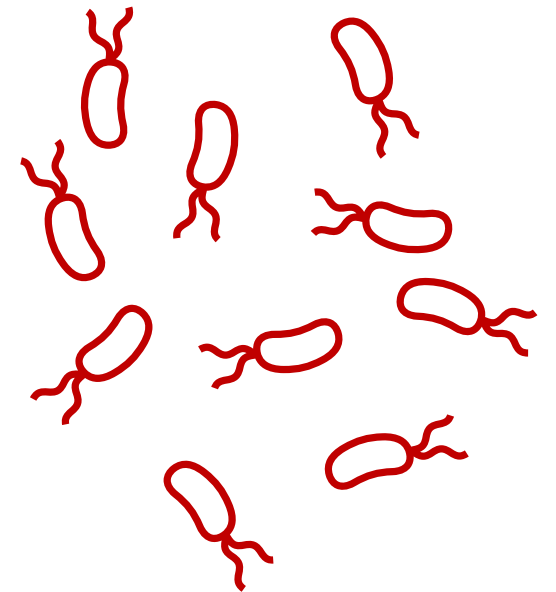
		 PC Positive Control	 NTC No Template Control	 Sample	 NEC Negative Extraction Control
Templates added to the real-time PCR	FAM Template		 Water	 DNA ready for PCR	 DNA extracted from sterile ATMP matrix
	ROX Template				
Read-outs	Target FAM Signal	 POSITIVE	 NEGATIVE	To be tested!	 NEGATIVE
	Internal Control ROX Signal	 Does not matter	 POSITIVE	To be tested!	 POSITIVE

 PCR grade Water  
 Positive Control  
 Internal Control

## Tips and tricks

- Extensive cleaning with chlorine-based agents
- Avoid cleaning with ethanol → Ethanol precipitates DNA
- Work carefully e.g. do not touch the lids of open tubes
- It is recommended to wear gloves and mask

Mycoplasma DNA  
is on our skin!



Simplifying Progress



Sartorius validation support

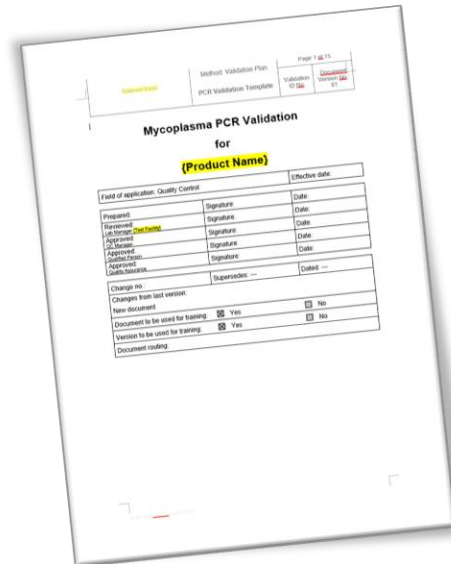
**SARTORIUS**

# Validation reports, templates & testing proposals



## ▪ Product Validation Report

- Microsart® ATMP Mycoplasma + Microsart® AMP Extraction



## ▪ Validation Template

- Microsart® ATMP Mycoplasma + Microsart® AMP Extraction



## ▪ Validation Proposal

- Standard matrix specific validation
- Individual support

## Further support for your validation

- Microsart® Validation Standard (10 CFU/Vial) & Microsart® Calibration Reagent (10<sup>8</sup> GC/Vial)
  - *Mycoplasma arginini*
  - *Mycoplasma orale*
  - *Mycoplasma gallisepticum*
  - *Mycoplasma pneumoniae*
  - *Mycoplasma synoviae*
  - *Mycoplasma fermentans*
  - *Mycoplasma hyorhinis*
  - *Acholeplasma laidlawii*
  - *Spiroplasma citri*
  - *Mycoplasma salivarium*

Non-viable CFU  
Standards!

DNA  
Standards!

Do you miss a species? Let us know!

# Application Notes

## Rapid, real-time PCR-based detection of microbial contaminations in high cell density Jurkat-, HPBMC-and CHO-cultures using Microsart® ATMP kits

In this study, we assessed the detection capability of Microsart® ATMP Extraction, combined with Microsart® ATMP Bacteria/Fungi/Mycoplasma assays, in high-density cell cultures.



Detection limits in different cell types

Cell Type Background	Microorganism Spike	Background Cells /mL (In 10 <sup>6</sup> )	Detection
CHO	99 CFU <i>B. subtilis</i>	19.0	Successful
CHO	10 CFU <i>M. arginini</i>	15.0 and 15.6 (two individual experiments)	Successful
CHO	10 CFU <i>M. orale</i>	15.5 and 16.3 (two individual experiments)	Successful
Jurkat	99 CFU <i>K. rhizophila</i>	10 to 40	Successful up to 25 x 10 <sup>6</sup> c/mL
Jurkat	50 CFU <i>C. albicans</i>	10 to 40	Successful up to 20 x 10 <sup>6</sup> c/mL
Jurkat	10 CFU <i>M. orale</i>	10 to 40	Not successful: PCR inhibition > 15 x 10 <sup>6</sup> c/mL No detection of Mycoplasma spike
Jurkat	10 CFU <i>M. synoviae</i>	10 to 35	Not successful: Partial PCR inhibition No detection of Mycoplasma spike
HPBMC	99 CFU <i>K. rhizophila</i>	10 to 40	Successful only up to 10 x 10 <sup>6</sup> c/mL
HPBMC	10 CFU <i>M. arginini</i>	15.0	Successful
HPBMC	10 CFU <i>M. orale</i>	19.1	Successful
HPBMC	99 CFU <i>P. aeruginosa</i>	20 and 25	Successful

Table 1: Results of Real-Time PCR detection of respective microbial spikes in varying cell types with respective cell densities.

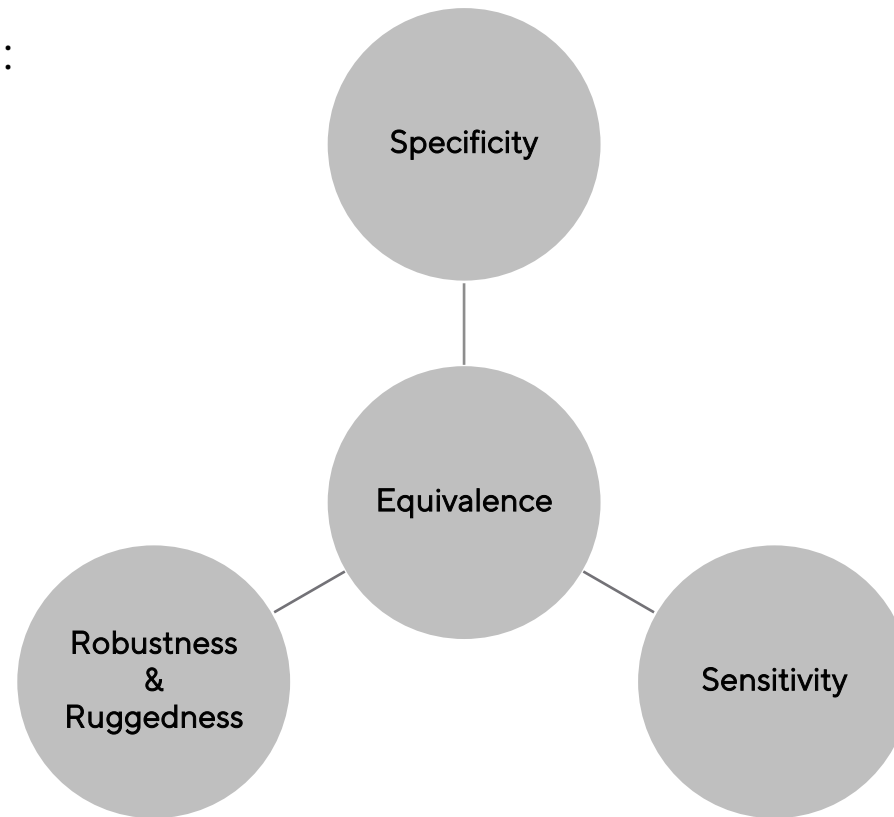
# Status quo regulatory landscape

Microbiological QC-Release testing			
Method	Mycoplasma	Bacteria	Fungi
Classical testing	USP<63>   EP 2.6.7 28 days	USP<71>   EP 2.6.1 Sterility testing 14 days	
real-time PCR-based	EP 2.6.7 (USP<1223>/EP 5.1.6)	USP<1071> EP 2.6.27 (USP<1223>/EP 5.1.6)	

# Validation overview

Regulatory guidance for validation:

- EP 2.6.7 (Mycoplasma)
- EP 2.6.21 (NAT)
- ICH Q2B
  
- USP<63>
- USP<1223>
- EP 5.1.6
- PDA TR 33



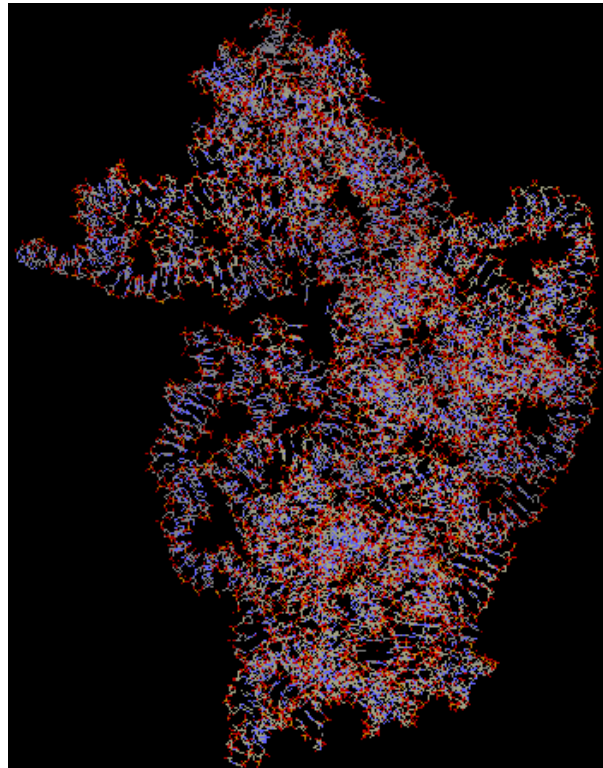
Including Guidance of the German Governmental Regulatory Agency (part of EMEA)

Paul-Ehrlich-Institut 



# *In silico* prediction by sequence alignment and blast

16S rDNA



<http://www.biochem.umd.edu/biochem/kahn/bchm465-01/ribosome/16SrRNA.html>

NIH U.S. National Library of Medicine

BLAST® » blastn suite

blastn blastp blastx tblastn tblastx

Enter Query Sequence

Enter accession number(s), gi(s), or FASTA sequence(s)

AGAGTTTGATCTGGCTCAG

Or, upload file

Job Title

Align two or more sequences

Choose Search Set

Database

Human genomic + transcript

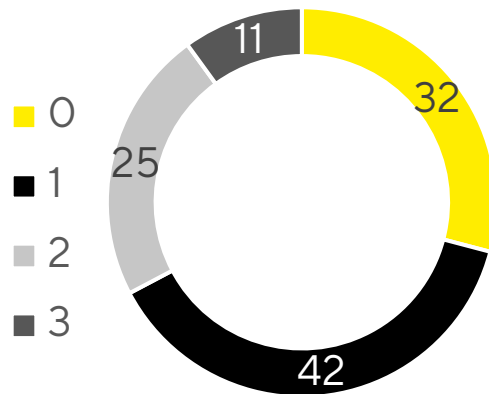
16S ribosomal RNA sequences (Bac

<https://www.ncbi.nlm.nih.gov/>



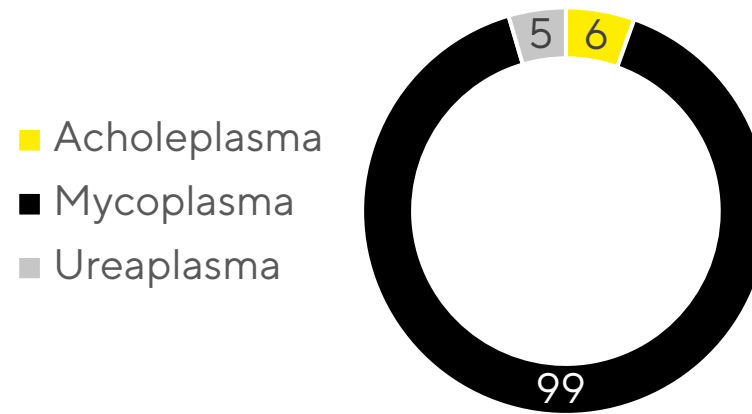
# Detection range

# Primer/Probe Mismatches

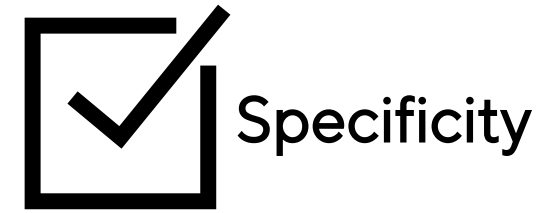


Acceptance criterion:  
≤ 3 nucleotides mismatch  
of primers and probe

Genus



At least **110 species** are detectable based on sequence alignment.



# Matrix effects

- Different matrices were tested for matrix effects and PCR inhibition
  - Typical cell culture media
  - Typical additives for cell culture media
  - Typical buffers

No.	Sample Matrix	Lot	Ct (FAM)	Ct (ROX)	Result
1	Hayflick Liquid Medium	107311	No Ct	35.01	negative
2	Frey Liquid Medium	107717	No Ct	32.39	negative
3	Dulbecco's MEM, bottle 1	1499W	No Ct	34.10	negative
4	Dulbecco's MEM, bottle 2	0407X	22.52	37.73	positive
5	RPMI Medium	0512X	No Ct	36.30	negative
6	PBS Dulbecco	0448W	No Ct	34.94	negative
7	anti-human-T-lymphocyte immune globulin, rabbit*	B 14 L-2	No Ct	34.24	negative
8	anti-human-T-lymphocyte immune globulin, rabbit*	B 13 L-1	No Ct	33.99	negative
9	FBS*	1231T	No Ct	35.13	negative
10	FBS*	11-3356	No Ct	34.65	negative
11	FBS Superior*	1201T	No Ct	34.61	negative
12	goat serum	B1060711	No Ct	33.96	negative
13	goat serum	A8070711	No Ct	33.78	negative
14	Tris buffer	118B1101	No Ct	33.67	negative
15	rabbit serum*	B03411-1243	No Ct	33.36	negative
16	donor horse serum	F11-2542	No Ct	34.11	negative
17	donor horse serum	0533W	No Ct	30.99	negative
18	calf serum*	11-3497	No Ct	32.46	negative
19	calf serum*	11-2875	No Ct	33.26	negative
20	horse serum*	10-2658	No Ct	34.32	negative
21	positive control	---	26.26	35.66	positive
22	NTC	---	No Ct	36.60	negative



16S rDNA sequencing revealed an *M. horvathii* contamination in Dulbecco's MEM bottle 2

No.	Sample Matrix	Ct (FAM)	Ct (ROX)	Result
1	Dulbecco's MEM, bottle 1, repeat 1	No Ct	32.76	negative
2	Dulbecco's MEM, bottle 1, repeat 2	No Ct	34.13	negative
3	Dulbecco's MEM, bottle 1, repeat 3	No Ct	34.28	negative
4	Dulbecco's MEM, bottle 1, repeat 4	No Ct	34.24	negative
5	Dulbecco's MEM, bottle 1, repeat 5	No Ct	33.90	negative
6	Dulbecco's MEM, bottle 2, repeat 1	No Ct	34.00	negative
7	Dulbecco's MEM, bottle 2, repeat 2	No Ct	34.70	negative
8	Dulbecco's MEM, bottle 2, repeat 3	No Ct	32.88	negative
9	Dulbecco's MEM, bottle 2, repeat 4	No Ct	33.43	negative
10	Dulbecco's MEM, bottle 2, repeat 5	No Ct	33.93	negative
11	positive control	28.47	34.37	positive
12	NTC	No Ct	35.34	negative
13	DNA extraction control	No Ct	36.40	negative

Second test on Dulbecco's MEM was contamination free and showed as well as the other matrices no matrix effects

For the tested matrices, no matrix effects were detected.

\* Due to confidentiality reasons no specifications can be provided for these matrices.



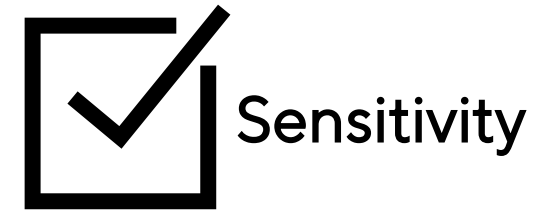
## Cross reactivity

- To exclude detection of other microorganisms or eukaryotic cells, the cross reactivity was assessed
- $\geq 0.1$  ng DNA/PCR for microorganisms and  $\geq 30$  ng for mammalian cells

Species	Results
<i>Clostridium acetobutylicum</i>	negative
<i>Lactobacillus acidophilus</i>	negative
<i>Streptococcus pneumoniae</i>	negative
Vero-B4	negative
Per.C6	negative
RK13	negative
CHO-K1	negative
Murine Genomic DNA	negative
Calf Thymus DNA	negative

For the tested organisms, **no cross reactivity** was detected.

# Limit of detection



EP 2.6.7

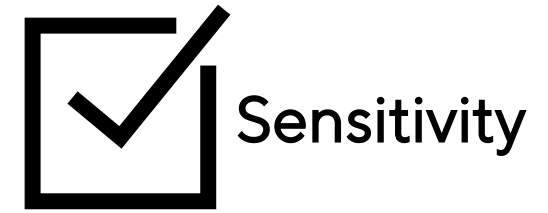
USP/EP required = 9



*Mycoplasma arginini*  
*Mycoplasma orale*  
*Mycoplasma gallisepticum*  
*Mycoplasma pneumoniae*  
*Mycoplasma synoviae*  
*Mycoplasma fermentans*  
*Mycoplasma hyorhinis*  
*Acholeplasma laidlawii*  
*Spiroplasma citri*

23/24 positive

Limit of detection (LOD<sub>95</sub>): 80, 40, 20, 10, 5, 2.5 CFU/ml



# Limit of detection




*Mycoplasma synoviae*

CFU/ml	80	40	20	10	5	2.5	NTC
	31.29	31.44	31.12	33.99	36.62	39.69	No Ct
	32.03	32.57	34.51	44.25	No Ct	33.78	
	31.94	32.02	33.28	39.00	33.39	35.09	
	30.14	33.07	30.80	33.77	35.26	39.36	No Ct
	30.69	31.24	30.85	33.72	No Ct	36.35	
	32.52	30.73	32.84	34.46	34.66	39.75	
	31.44	33.24	32.94	33.95	No Ct	36.92	No Ct
	31.27	32.38	36.52	33.12	38.97	36.72	
	29.34	33.88	34.26	No Ct	33.90	36.86	
	32.00	33.32	32.11	39.01	33.64	32.85	No Ct
	31.43	31.98	34.76	30.82	34.94	37.47	
	33.18	31.48	35.30	37.10	30.73	31.42	
	32.20	33.53	33.45	36.00	34.03	39.85	No Ct
	32.07	33.17	34.68	37.45	34.26	37.35	
	32.26	31.91	34.49	36.09	41.95	No Ct	
	29.82	32.51	31.45	33.07	39.94	33.94	No Ct
	32.54	33.48	31.96	33.44	31.55	32.08	
	30.38	33.88	32.56	32.98	34.10	33.19	
	32.92	34.47	34.45	36.28	32.61	36.36	No Ct
	33.86	36.67	37.23	36.72	35.16	38.16	
	33.90	33.95	35.18	35.89	34.98	39.63	
	31.63	32.52	32.16	31.06	33.58	29.66	No Ct
	30.75	31.36	32.00	29.83	33.52	33.00	
	33.32	28.08	32.42	28.90	34.03	32.52	
Mw	31.79	32.62	33.39	34.82	34.85	35.74	
STABWN	1.184	1.574	1.716	3.305	2.580	2.963	
positive	24	24	24	23	20	23	
total	24	24	24	24	24	24	

Species (CFU-based)	Acceptance criterion	LOD <sub>95</sub> (CFU/mL)
<i>Mycoplasma arginini</i>	23/24	10
<i>Mycoplasma orale</i>	23/24	10
<i>Mycoplasma gallisepticum</i>	23/24	10
<i>Mycoplasma pneumoniae</i>	23/24	≤ 5
<i>Mycoplasma synoviae</i>	23/24	10
<i>Mycoplasma fermentans</i>	23/24	10
<i>Mycoplasma hyorhinitis</i>	23/24	≤ 5
<i>Acholeplasma laidlawii</i>	23/24	10
<i>Spiroplasma citri</i>	23/24	≤ 5



# Device comparability

Cycler	Microsart® ATMP Mycoplasma
 Thermo Fischer ABI Prism 7500	<b>PASSED</b>
 Agilent Mx3005p	<b>PASSED</b>
 Qiagen Rotor-Gene 6000	<b>PASSED</b>

... by today many more devices are used by our customers.

# Validation overview



- Sensitivity
  - LOD<sub>95</sub> – limit of detection
  - Culture media comparison
- Specificity
  - Sequence alignment
  - Sample matrix effects
  - Cross reactivity
  - Mollicutes detection range
  - Identification of unspecific amplification
- Robustness
  - Cell culture material
  - Real-time PCR cycler compatibility



Simplifying Progress



Customer Validation Data  
Service Lab Labor Dr. Quade



SARTORIUS



# Microsart® ATMP Mycoplasma Real-time PCR – Matrix specific Validation with automated DNA extraction

- Service Lab Labor Dr. Quade provides mycoplasma testing service for a cell therapy manufacturer (release testing of a T-cell-based product)
  - Samples
    - transduced **human T lymphocytes** at a concentration of **12.5 million cells per mL** in cultivation medium
  - Mycoplasma contamination detection system
    - automated DNA extraction using MagnaPure Compact Instrument
    - Microsart® ATMP Mycoplasma
  - Spikes
    - Sartorius Microsart® Validation Standards (10 CFU)
    - included species: *M. orale*, *M. fermentans*, *M. pneumoniae*, *A. laidlawii*
  - Acceptance criterion
    - 23 positive signals out of 24 replicates (each species)



# Microsart® ATMP Mycoplasma Real-time PCR – Matrix specific Validation with automated DNA extraction

Table 1: *Mycoplasma pneumoniae*

Sample spiked with 10 CFU/mL	Result CP 19.11.2019	Result CP 20.11.2019	Result CP 21.11.2019	Inter-Assay CV%
Sample 1	29.93	29,23	32,99	6,51
Sample 2	29.85	29,60	32,85	5,88
Sample 3	29,33	29,27	32,27	5,66
Sample 4	30.00	29,63	32,47	5,03
Sample 5	29.87	29,58	32,24	4,77
Sample 6	30.08	29,50	32,45	5,09
Sample 7	29.73	29,79	32,09	4,41
Sample 8	29,87	30,03	31,92	3,73
Intra-Assay CV%	0.77	0,88	1.12	

CP: Crossing Point, Coefficient of Variance in %, Negative samples showed CP value >40 or a plane straight line



# Microsart® ATMP Mycoplasma Real-time PCR – Matrix specific Validation with automated DNA extraction

Table 2: *Mycoplasma fermentans*

Sample spiked with 10 CFU/mL	Result CP 20.11.2019	Result CP 21.11.2019	Result CP 22.11.2019	Inter-Assay CV%
Sample 1	33,60	33,15	31,29	3,75
Sample 2	33,65	32,67	31,57	3,19
Sample 3	33,35	32,83	31,50	2,93
Sample 4	32,91	33,22	31,61	2,62
Sample 5	33,44	32,94	31,48	3,12
Sample 6	32,99	32,79	31,55	2,40
Sample 7	32,80	33,12	31,80	2,11
Sample 8	33,16	30,00	31,56	5,00
Intra-Assay CV%	0,97	3,27	0,45	

CP: Crossing Point, Coefficient of Variance in %, Negative samples showed CP value >40 or a plane straight line



# Microsart® ATMP Mycoplasma Real-time PCR – Matrix specific Validation with automated DNA extraction

Table 3: *Mycoplasma orale*

Sample spiked with 10 CFU/mL	Result CP 20.11.2019	Result CP 21.11.2019	Result CP 22.11.2019	Inter-Assay CV%
Sample 1	31,53	32,47	31,64	1,61
Sample 2	31,56	32,68	31,93	1,78
Sample 3	31,68	33,02	31,21	2,94
Sample 4	31,21	32,36	31,84	1,81
Sample 5	31,19	32,48	31,27	2,28
Sample 6	31,20	32,86	31,54	2,75
Sample 7	31,23	32,14	31,43	1,51
Sample 8	31,00	32,47	32,04	3,50
Intra-Assay CV%	0.75	1.06	0.97	

CP: Crossing Point, Coefficient of Variance in %, Negative samples showed CP value >40 or a plane straight line



# Microsart® ATMP Mycoplasma Real-time PCR – Matrix specific Validation with automated DNA extraction

Table 4: *Acholeplasma laidlawii*

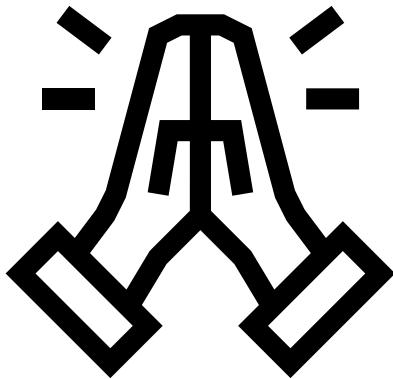
Sample spiked with 10 CFU/mL	Result CP 20.11.2019	Result CP 21.11.2019	Result CP 22.11.2019	Inter-Assay CV%
Sample 1	31,46	28,61	31,06	5,08
Sample 2	31,28	28,17	30,73	5,52
Sample 3	31,07	29,00	31,27	4,13
Sample 4	31,52	29,00	30,93	4,32
Sample 5	31,12	31,29	31,01	0,45
Sample 6	30,97	31,33	30,94	0,70
Sample 7	31,38	31,55	31,15	0,64
Sample 8	30,74	31,44	30,98	1,15
Intra-Assay CV%	0.85	4.9	0.52	

CP: Crossing Point, Coefficient of Variance in %, Negative samples showed CP value >40 or a plane straight line



# Microsart® ATMP Mycoplasma Real-time PCR – Matrix specific Validation with automated DNA extraction

- Matrix specific validation was **successful**
- The transduced **human T lymphocytes** at a concentration of **12.5 million cells per mL** in cultivation medium are suitable for release testing using automated extraction and the Microsart® ATMP Mycoplasma kit



# SARTORIUS

## Simplifying Progress

Microsart® AMP Extraction  
Microsart® AMP Mycoplasma

2022





# Mycoplasma contamination detection

- Real-time PCR allows detection of Mycoplasma
  - In 3 h
  - Down to 5-10 CFU/ml
- Validated combination
  - In accordance with EP 2.6.7 and USP 63
- Support
  - *Product Validation Report* containing all experimental details
  - *Matrix Validation Template* containing detailed information for the customer specific matrix validation
  - Technical support during matrix validation process



Microsart® AMP Extraction



Microsart® AMP Mycoplasma

# Workflow Mycoplasma contamination detection

- Concentration of the matrix to be tested
  - Vivaspin® 6 or Vivaspin® 20
  - Volume is reduced to 200 µl sample
- DNA isolation using the column-based Microsart® AMP Extraction kit
  - Columns allow to isolate DNA from the whole 200 µl ATMP sample (supernatant & cells)
- Real-time PCR using the Microsart® AMP Mycoplasma kit
  - PCR contains 50 µl sample DNA → complete analysis of 20 ml sample
  - Taq-Man® System → reduce false-positive signals
  - Duplex assay → reduce false-negative signals
  - Universal assay for different real-time PCR cycler → High volume thermo cycling, FAM™ and ROX™
  - High stability & no freezing → Lyophilized reagents



20 mL



6 mL

100.000 MWCO



# RESEARCH



# SARTORIUS

## Simplifying Progress

### Microsart® RESEARCH Mycoplasma

2022



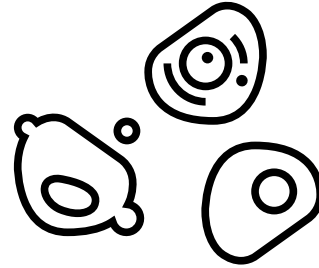
# In-process contamination control of Mycoplasma

- Key advantages

- Very robust towards inhibitors
- No prior DNA extraction required
- Internal control DNA included in real-time PCR master mix
- One step preparation

→ Quick 'n' Dirty for process monitoring

- Taq-Man® System → reduce false-positive signals
- Duplex assay → reduce false-negative signals
- Universal assay for different real-time PCR cycler → FAM™ and ROX™
- High stability & no freezing → Lyophilized reagents



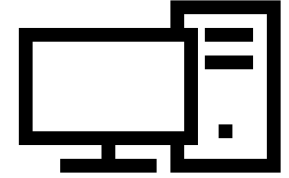
ATMP  
(CHO, HEK,...)



2 µl



23 µl  
RESEARCH  
master mix



Detection in  
real-time PCR cycler:  
Contamination? Yes/No



Microsart® RESEARCH  
Mycoplasma

# Thank you.



PCR@Sartorius.com

**SARTORIUS**