



Jena Bioscience, with over 25 years of experience, is a leading provider of inovative and high-quality reagents and customized services in the life science field.



**Bulk – Tailored to your scale.**

Ready-made or customized life science reagents (liquids, lyo-ready or lyophilized) in large quantities, with personalized packaging and scientific support.



**OEM – Your products, your way.**

Exclusivity with customized formulations, packaging and branding under strict confidentiality.

**Contact our Molecular Biology Experts:**

Phone: +49(0)3641 - 6285 000  
Mail: molbio@jenabioscience.com



**Dr. Bürk Schäfer**  
Head of Molecular Biology Dept.



**Dr. Juliane Buschmann**  
Business Development Manager



## LIQUID MIXES FOR LYOPHILIZATION

Optimized for lyophilization in your own laboratory

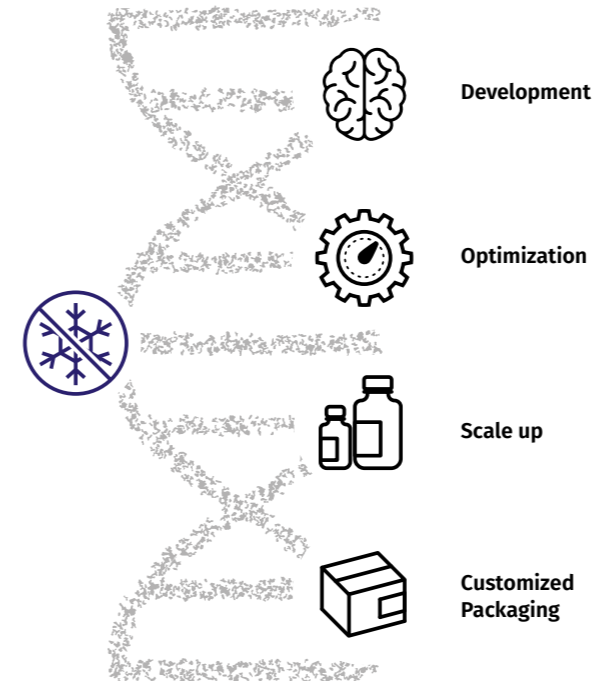


Cat.-No.	Amount	Conc.	Reactions
<b>qPCR SybrMaster Lyophilisate</b>			
PCR-189-1ML	1 ml	2.5 ×	125 reactions × 20 µl
PCR-189-10ML	10 ml	2.5 ×	1,250 reactions × 20 µl
<b>qPCR ProbesMaster Lyophilisate</b>			
PCR-188-1ML	1 ml	2.5 ×	125 reactions × 20 µl
PCR-188-10ML	10 ml	2.5 ×	1,250 reactions × 20 µl
<b>Liquid Direct qPCR SybrMaster Lyophilisate</b>			
PCR-187-1ML	1 ml	2.5 ×	125 reactions × 20 µl
PCR-187-10ML	10 ml	2.5 ×	1,250 reactions × 20 µl
<b>Liquid Direct qPCR ProbesMaster Lyophilisate</b>			
PCR-186-1ML	1 ml	2.5 ×	125 reactions × 20 µl
PCR-186-10ML	10 ml	2.5 ×	1,250 reactions × 20 µl

## CUSTOMIZATION

Looking for freeze-dried mixes beyond our standard repertoire?

We are your reliable partner with all technical advice and know how for developing lyophilisates according to your requirements.



Contact us for customized lyophilization service:  
[molbio@jenabioscience.com](mailto:molbio@jenabioscience.com)



# LYOPHILIZED REAGENTS FROM JENA BIOSCIENCE

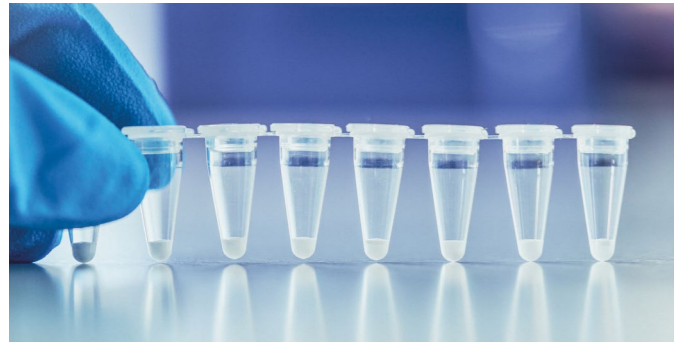


Ready-to-Use – Pre-aliquoted assays  
No fridge required – Lowest risk of contamination – Most simple assay set-up



## qPCR

All our lyophilized master mixes contain polymerases and nucleotides of highest quality and are optimized for maximum sensitivity as well as specificity



Lyophilisates based on the detection with **SYBR® Green** are perfect for setting up singleplex reactions without the need of extensive optimization.

For **TaqMan® Detection** including duplex or multiplex assay setups, ProbesMaster Lyophilisates are the first choice.

Cat.-No.	Amount
<b>qPCR SybrMaster Lyophilisate</b>	
PCR-173S	192 reactions × 20 µl
PCR-173L	960 reactions × 20 µl
<b>qPCR ProbesMaster Lyophilisate</b>	
PCR-156S	192 reactions × 20 µl
PCR-156L	960 reactions × 20 µl

## DIRECT qPCR

Save on time- and money-consuming DNA extraction – perfect for Point-of-Care applications

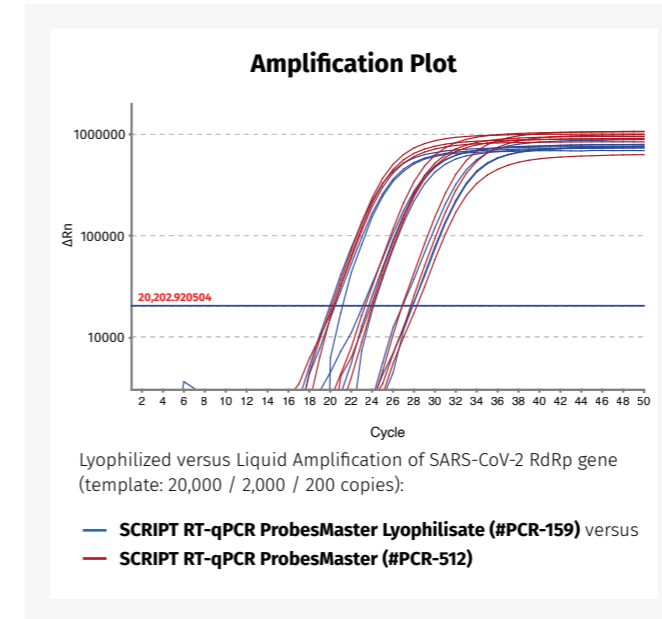


- Automatable for high throughput
- Reduce DNA preparation time by 70–90 %
- Time & cost efficient
- Minimize sample loss
- Less plastic consumables

Cat.-No.	Amount
<b>Direct qPCR SybrMaster Lyophilisate</b>	
PCR-177S	192 reactions × 20 µl
PCR-177L	960 reactions × 20 µl
<b>Direct qPCR ProbesMaster Lyophilisate</b>	
PCR-176S	192 reactions × 20 µl
PCR-176L	960 reactions × 20 µl

## RT-qPCR

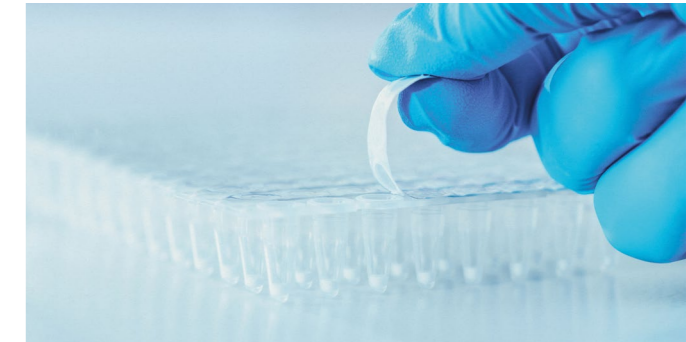
Ready-to-use lyophilisates come as complete master mixes in a dry, room temperature stable format



Cat.-No.	Amount
<b>SCRIPT RT-qPCR SybrMaster Lyophilisate</b>	
PCR-169S	192 reactions × 20 µl
PCR-169L	960 reactions × 20 µl
<b>SCRIPT RT-qPCR ProbesMaster Lyophilisate</b>	
PCR-159S	192 reactions × 20 µl
PCR-159L	960 reactions × 20 µl

## LAMP

Isothermal Amplification – Revolutionize the way for detecting DNA



Fast, simple and cost-effective: generate results without the need of expensive instruments.

**LAMP assays for fast (10–20 min) and ultra-fast (5–10 min) detection of DNA**

Cat.-No.	Amount
<b>Saphir Bst GreenMaster Lyophilisate</b>	
PCR-398S	192 reactions × 20 µl
PCR-398L	960 reactions × 20 µl
<b>Saphir Bst Turbo GreenMaster Lyophilisate</b>	
PCR-395S	192 reactions × 20 µl
PCR-395L	960 reactions × 20 µl