



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



Your partner for freeze-drying

LIQUID MIXES FOR LYOPHILIZATION

Optimized for lyophilization in your own laboratory

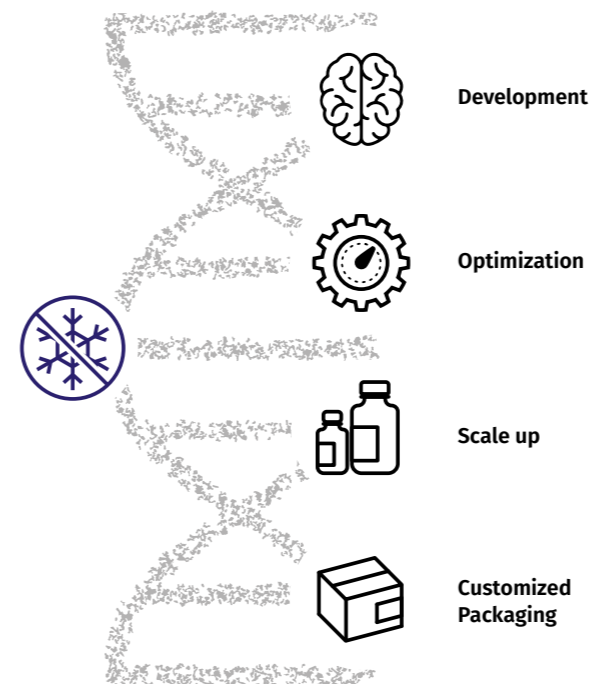


Cat.-No.	Amount	Conc.	Reactions
Liquid qPCR GreenMaster Lyophilisate			
PCR-189-1ML	1 ml	2.5 ×	125 reactions × 20 µl
PCR-189-10ML	10 ml	2.5 ×	1,250 reactions × 20 µl
Liquid qPCR ProbesMaster Lyophilisate			
PCR-188-1ML	1 ml	2.5 ×	125 reactions × 20 µl
PCR-188-10ML	10 ml	2.5 ×	1,250 reactions × 20 µl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



LYOPHILIZED REAGENTS FROM JENA BIOSCIENCE

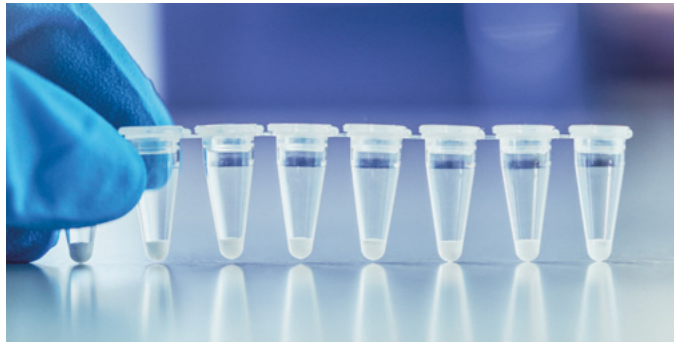


Ready-to-Use – Pre-aliquoted assays
No fridge required – Lowest risk of contamination – Most simple assay set-up



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 a **green-fluorescent dye** 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
qPCR a GreenMaster dye Lyophilisate	
PCR-173S	192 reactions × 20 µl
PCR-173L	960 reactions × 20 µl
qPCR ProbesMaster Lyophilisate	
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

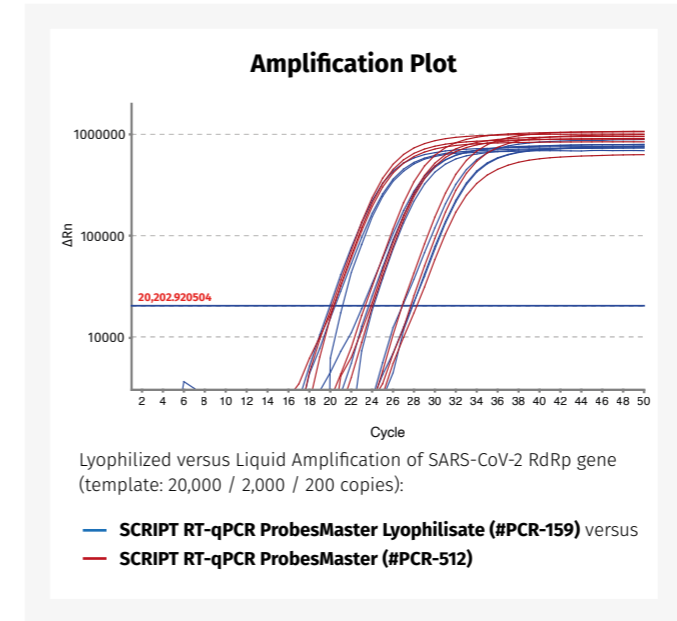


- Automatable for high throughput
- Reduce DNA preparation time by 70–90 %
- Time & cost efficient
- Minimize sample loss
- Less plastic consumables

Cat.-No.	Amount
Direct qPCR GreenMaster Lyophilisate	
PCR-177S	192 reactions × 20 µl
PCR-177L	960 reactions × 20 µl
Direct qPCR ProbesMaster Lyophilisate	
PCR-176S	192 reactions × 20 µl
PCR-176L	960 reactions × 20 µl

RT-qPCR

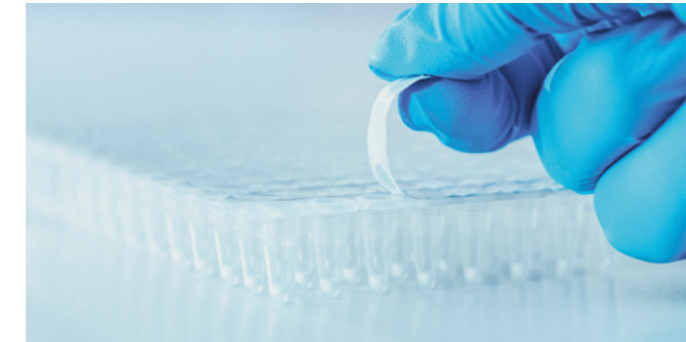
Ready-to-use lyophilisates come as complete master mixes in a dry, room temperature stable format



Cat.-No.	Amount
SCRIPT RT-qPCR GreenMaster Lyophilisate	
PCR-169S	192 reactions × 20 µl
PCR-169L	960 reactions × 20 µl
SCRIPT RT-qPCR ProbesMaster Lyophilisate	
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
Saphir LAMP GreenMaster Lyophilisate	
PCR-398S	192 reactions × 20 µl
PCR-398L	960 reactions × 20 µl
Saphir LAMP Turbo GreenMaster Lyophilisate	
PCR-395S	192 reactions × 20 µl
PCR-395L	960 reactions × 20 µl