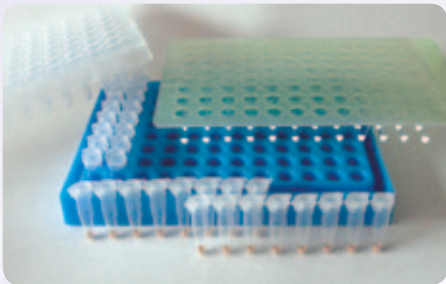


PCR Lyophilisates

PCR Lyophilisates – Preloaded and Stable at Room Temperature



Jena Bioscience lyophilized mastermixes

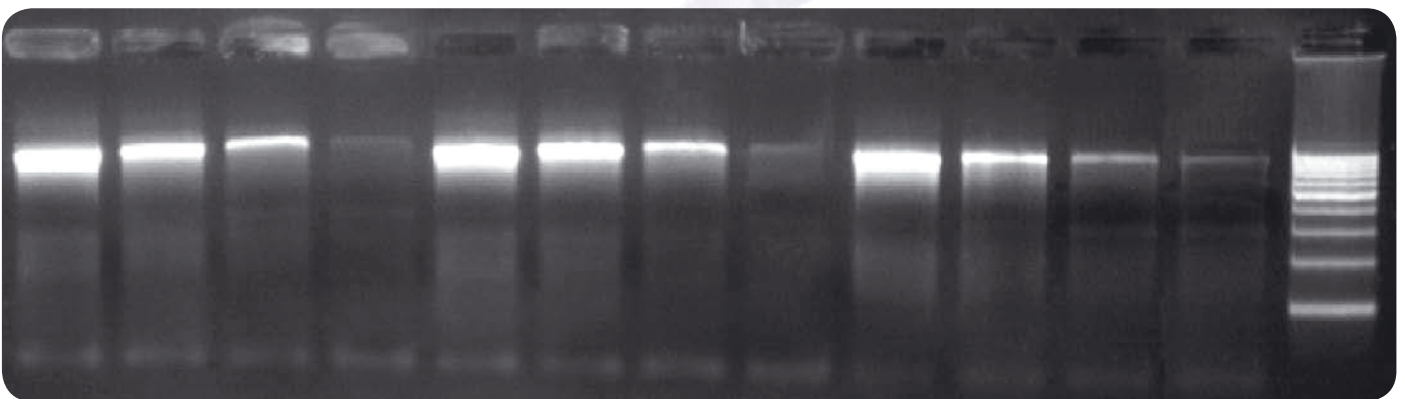
- are available for standard PCR, real-time-PCR and reverse transcription
- provide “point of care” suitability due to stability at 30°C for at least 6 months
- combine highest performance with convenience of use
- guarantee highest DNA yields
- minimize risk of pipetting errors or contamination

Ready-to-Use Lyophilisates are delivered in PCR reaction tube strips and as 96-well plates preloaded with a complete master mix in a dry, room temperature stable format. There is no need for freezing, thawing or pipetting on ice. Simply fill up the vials with template DNA, primer and PCR-grade water and run the PCR in any thermocycler.

Provide us your primer sequence and we deliver preloaded PCR tubes and plates ready-made for your special application.

Upon request we prepare lyophilisates comprising primers, polymerase, dNTPs and reaction buffer. All you still have to do is adding template DNA, filling up with water and starting the cycler!

Taq Master Lyophilisates show full functionality after storage at 30°C for 6 months



100 pg	50 pg	25 pg	10 pg	100 pg	50 pg	25 pg	10 pg	100 pg	50 pg	25 pg	10 pg	500 bp Ladder
Storage at 20°C for 6 months				Storage at 30°C for 6 months				Storage at -20°C for 6 months				

Assay: Amplification of a 4 kb fragment from lambda DNA, template dilution series • Cycling: 2 min 95°C, (10 sec 95°C, 20 sec 60°C, 4 min 72 °C) x 30, 4 min 72°C



Jena Bioscience GmbH
Loebstedter Str. 80
07749 Jena
Germany

Phone +49 (0)3641-62 85 000
Fax +49 (0)3641-62 85 100
www.jenabioscience.com
info@jenabioscience.com

 Jena Bioscience
www.jenabioscience.com

PCR Lyophilisates

Ready-to-Use Lyophilisates

Product	Cat.-No.	Size	Price
qPCR ProbesMaster Lyophilisate	PCR-156S-8TS	12 strips (96 reactions)	150,- Euro
	PCR-156L-8TS	60 strips (480 reactions)	600,- Euro
qPCR GreenMaster Lyophilisate	PCR-157S-8TS	12 strips (96 reactions)	150,- Euro
	PCR-157L-8TS	60 strips (480 reactions)	600,- Euro
Red Load Taq Master Lyophilisate	PCR-151S-8TS	12 strips (96 reactions)	90,- Euro
	PCR-151L-8TS	60 strips (480 reactions)	360,- Euro
Red Load Taq Master Lyophilisate / high yield	PCR-154S-8TS	12 strips (96 reactions)	90,- Euro
	PCR-154L-8TS	60 strips (480 reactions)	360,- Euro
Taq Master Lyophilisate	PCR-152S-8TS	12 strips (96 reactions)	90,- Euro
	PCR-152L-8TS	60 strips (480 reactions)	360,- Euro
Taq Master Lyophilisate / high yield	PCR-155S-8TS	12 strips (96 reactions)	90,- Euro
	PCR-155L-8TS	60 strips (480 reactions)	360,- Euro
Hot Start Master Lyophilisate	PCR-153S-8TS	12 strips (96 reactions)	120,- Euro
	PCR-153L-8TS	60 strips (480 reactions)	480,- Euro
One-Step RT-PCR Master Lyophilisate	PCR-159S-8TS	12 strips (96 reactions)	420,- Euro
	PCR-159L-8TS	60 strips (480 reactions)	1.680,- Euro
RT Master Lyophilisate	PCR-158S-8TS	12 strips (96 reactions)	280,- Euro
	PCR-158L-8TS	60 strips (480 reactions)	1.120,- Euro

Please contact us for inquiries and further information or request your customized quotation.

Find out more details at our website
www.jenabioscience.com



Jena Bioscience GmbH
 Loebstedter Str. 80
 07749 Jena
 Germany

Phone +49 (0)3641-62 85 000
 Fax +49 (0)3641-62 85 100
www.jenabioscience.com
info@jenabioscience.com



Jena Bioscience
www.jenabioscience.com