



**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



**SINGLE COMPONENTS**

**Polymerases, nucleotides, buffers and additives in individual quantities or fillings**

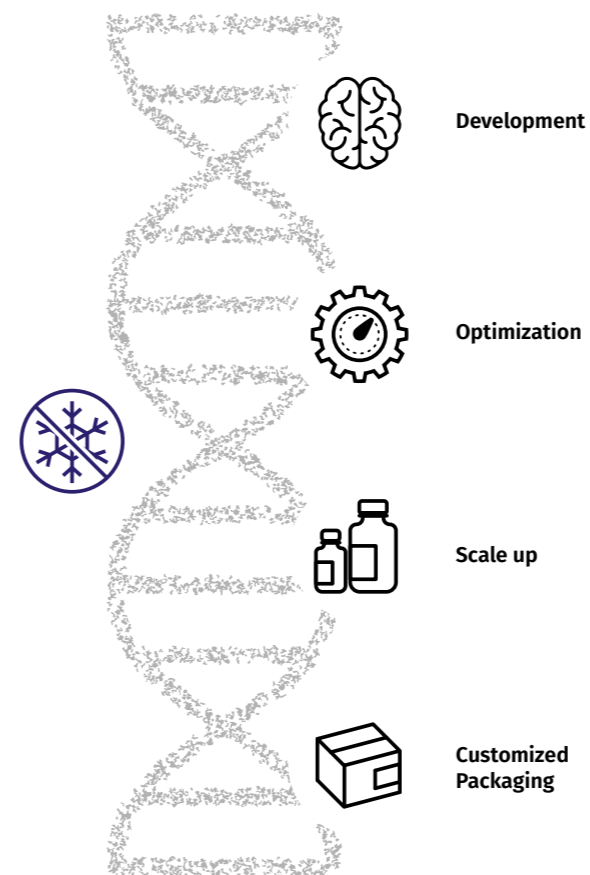
- Polymerases & Enzymes**
  - DNA & RNA polymerases
  - Thermolabile UNG
  - RNase inhibitor
- Lyo-Ready Components**
  - Glycerol-free polymerases & enzymes
  - Highly concentrated reagents for lyophilization
- Nucleotides**
  - In-house manufactured
  - Ultrapure (> 99.0 %)
  - Singles, bundles or mixtures
- Buffer, Water & Supplements**
  - Pre-mixed buffers
  - PCR-grade water
  - Gel loading buffer
  - DNA staining

- Complete control over your product
- Product optimization
- Change in formulation
- Kit configuration

Your Partner in  
Molecular Biology

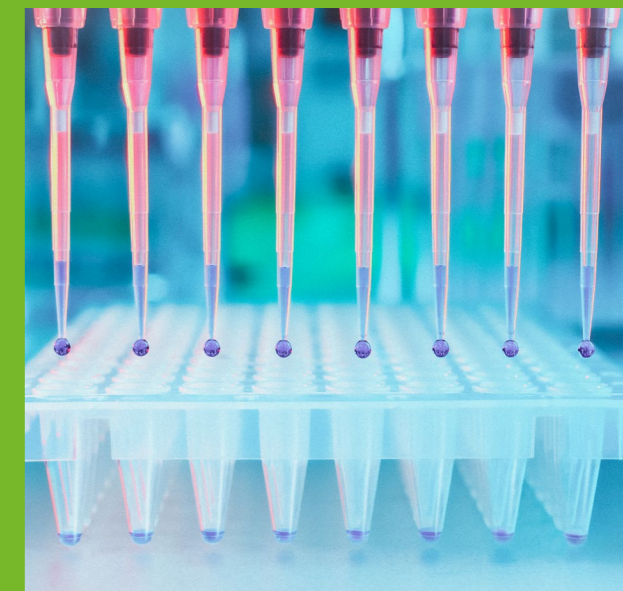
**LYOPHILIZATION**

Lyophilisates according to your requirements



Contact us for customized lyophilization service:  
**molbio@jenabioscience.com**

**YOUR PARTNER  
IN MOLECULAR  
BIOLOGY**



**Ready-to-Use & Customized Kits  
Liquid & Lyophilized – OEM Solutions  
Lab to Bulk Scale**



## DNA/RNA PREPARATION & CLEANUP

### DNA Extraction Kits

Efficient extraction kits based on columns or solutions, suitable for various sources

### RNA Isolation Kits

Rapid and effective isolation of highly purified RNA

### Cleanup Kits

Remove impurities, contaminants and enzymatic inhibitors

### Plasmid Preps

Column based isolation of plasmid or cosmid DNA

### Genomic DNA, plasmid DNA, total RNA or viral RNA from:



Tissue



Swabs



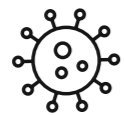
Blood



Bacteria



Yeast



Viruses

## CLASSICAL PCR

Product range for End point PCR

- High quality thermophilic polymerases
- Ultrapure nucleotides

### Master Mixes

- End point PCR
- Multiplex
- Direct gel loading
- Hot Start

### Polymerases & Core Kits

- Taq, Hot Start, High Fidelity, Phusion
- Proof-reading
- Aptamer inhibited
- Antibody blocked

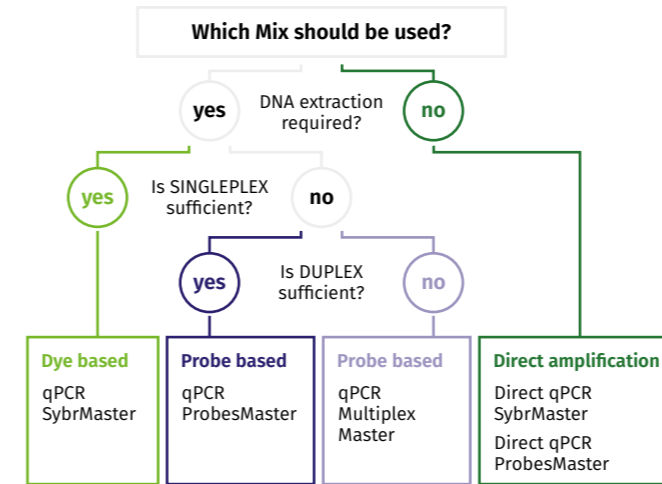
### DNA Ladders

- Log scale
- Linear scale
- Fluorescent
- Digests of DNA

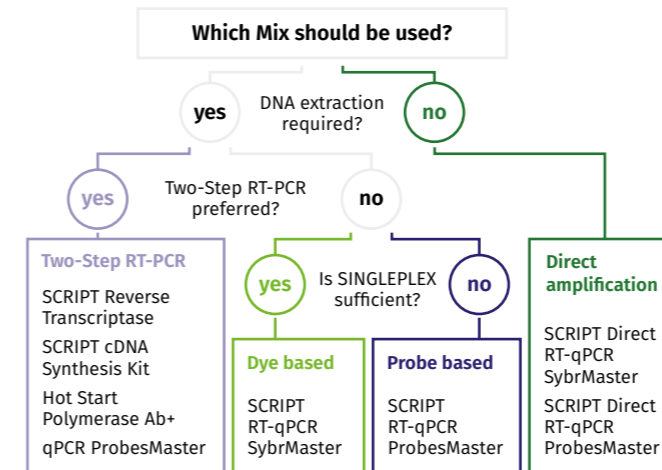
### Chromatography

- Agarose
- Loading buffer
- DNA staining

## qPCR



## RT-qPCR



## DIRECT AMPLIFICATION

No need for time-consuming DNA extraction

- Cut down total time for DNA analysis
- Automatable for high throughput
- Perfect for point-of-care applications
- Time & cost efficient
- Less plastic waste

## ISOTHERMAL AMPLIFICATION & LAMP

Loop mediated isothermal amplification

**Isothermal amplification**, with the standout technique known as **LAMP** (Loop-mediated Isothermal Amplification) is revolutionizing the way we analyze DNA and RNA.

LAMP is performed at a constant temperature (60–65°C), so no thermal cycling is required. Optimized Bst enzymes and master mixes generate an amplification factor of up to 10<sup>9</sup>, comparable to 30 cycles in a PCR assay. This allows the detection of a target gene within 10–30 minutes.

- Saphir Bst Polymerase
- Saphir LAMP GreenMaster
- Saphir LAMP GreenMaster Lyophilisate