

# OUR BENEFITS FOR SCIENTISTS FROM PARADISE

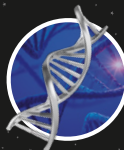
## CLICK CHEMISTRY

Convenient bioorthogonal molecule coupling (e.g. DNA & fluorescent dye)



## MOLECULAR BIOLOGY

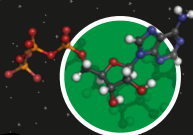
- PCR, qPCR & RT-PCR
- Cloning & mutagenesis
- DNA/RNA preparation & cleanup



## NUCLEOTIDES & NUCLEOSIDES

Natural & modified

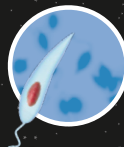
- NTPs
- dNTPs
- ddNTPs
- Nucleosides



## LEXSY EXPRESSION

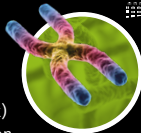
Eukaryotic protein expression solutions

- *In vivo* and *in vitro* system
- Custom protein expression



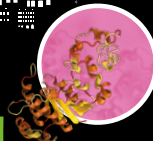
## PROBES & EPIGENETICS

- Non-radioactive biomolecule (DNA, RNA, protein) labeling (fluorescent, Hapten, Click)
- DNA & RNA (De)Methylation
- Protein (De)Acetylation



## PROTEINS

> 500 functionally active and highly purified recombinant & native proteins



## Jena only!

### Your BuyLocal Benefits:



7,5% discount on all products  
(except Oligos)



Fast delivery: we're just around  
the corner!



Same day pickup for all in-stock  
items



Individual mixes, concentrations  
or packaging



Eco-friendly due to short trans-  
portation routes: we are certified  
for our environmental manage-  
ment (EMAS certificate)



Easy ordering via webshop

## Any questions?



**Dr. Barbara Zschoernig**

Click Chemistry,  
Probes & Epigenetics

**Dr. Andreas Licht**

LEXSY Expression

**Christin Reuter**

Crystallography  
& Cryo-EM

**Dr. Buerk Schaefer**

Molecular Biology  
& Proteins

**Dr. Thomas Waldbach**

Nucleotides &  
Nucleosides

Get in touch with our product development scientists.



**Jena Bioscience**  
Building Blocks of Life

**Jena Bioscience GmbH**  
Loebstedter Str. 71  
07749 Jena, Germany

Phone +49(0)3641-6285 000  
Fax +49(0)3641-6285 100  
buylocal@jenabioscience.com  
[www.jenabioscience.com](http://www.jenabioscience.com)