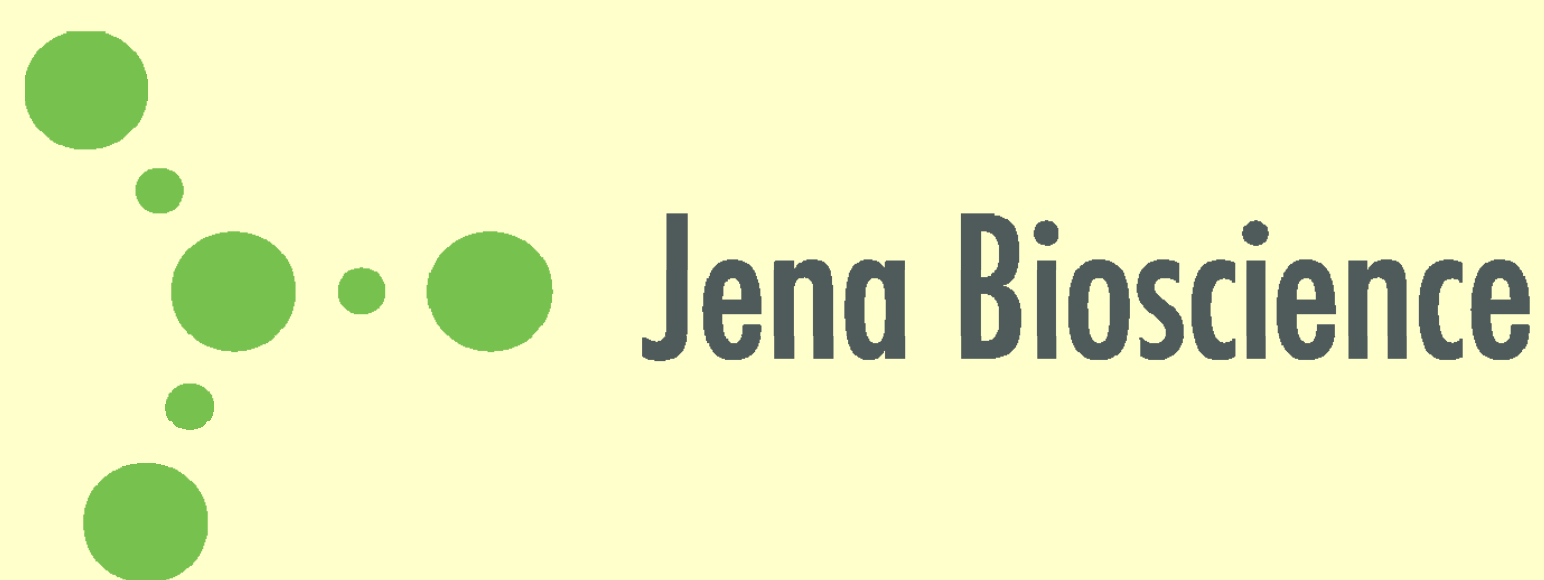


# The LEXSY Technology - a valuable tool in *Kinetoplastida*

R. Breitling<sup>1</sup>, S. Kushnir<sup>2</sup>, Y. Heinekamp<sup>1</sup>, I. Cirstea<sup>2</sup>, S. Mureev<sup>2,3</sup>, O. Kovtun<sup>2,3</sup> and K. Alexandrov<sup>2,3</sup>

<sup>1</sup> Jena Bioscience GmbH, Jena, Germany; <sup>2</sup> Max-Planck-Institut für molekulare Physiologie, Dortmund, Germany

<sup>3</sup> Institute for Molecular Biosciences, University of Queensland, Australia



## ABSTRACT

We have engineered the non-pathogenic parasite of lizard *Leishmania tarentolae* into a flexible and efficient protein expression platform called LEXSY. This biosafety 1 host system can be used for expression of target proteins of other *kinetoplastida*, making use of the close relationship of cellular machineries including protein folding and modification.

On the other hand the LEXSY vectors are functionally in other *Leishmania* species and were used successfully in *L. amazonensis*, *L. donovani*, *L. infantum*, *L. major*, *L. mexicana* and also in the plant parasite *Phytomonas serpens*.

Two principal LEXSY architectures yield constitutive or inducible expression. The constitutive version is based on integration of expression cassettes into the 18S rRNA gene (*ssu*) for transcription by the strong RNA polymerase I of the host.

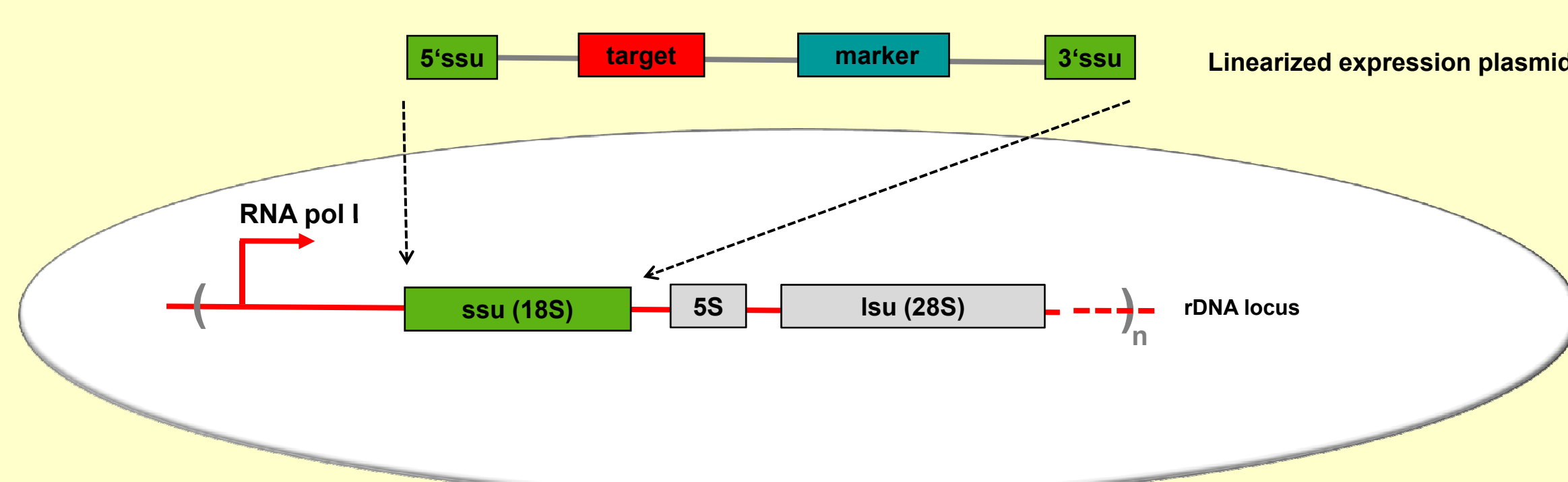
Four alternative antibiotic selection markers permit integration of up to four expression cassettes into the *ssu* locus, useful e.g. for expression of multisubunit proteins.

The inducible version employs T7 RNA Polymerase and TET repressor, expressed from the *ssu* locus of an engineered host. Target gene cassettes under control of a T7 promoter TET operator assembly can be integrated into  $\beta$ -tubulin ( *$\beta$ -tub*) or ornithine decarboxylase (*odc*) loci. Alternatively, expression cassettes can be positioned on circular or linear episomes with telomeric ends. The use of a fusion of the *bleo* resistance and *cherry* fluorescence genes provides a tool for selection and screening.

The LEXSY toolbox has recently been complemented by a cell-free version for expression of recombinant proteins.

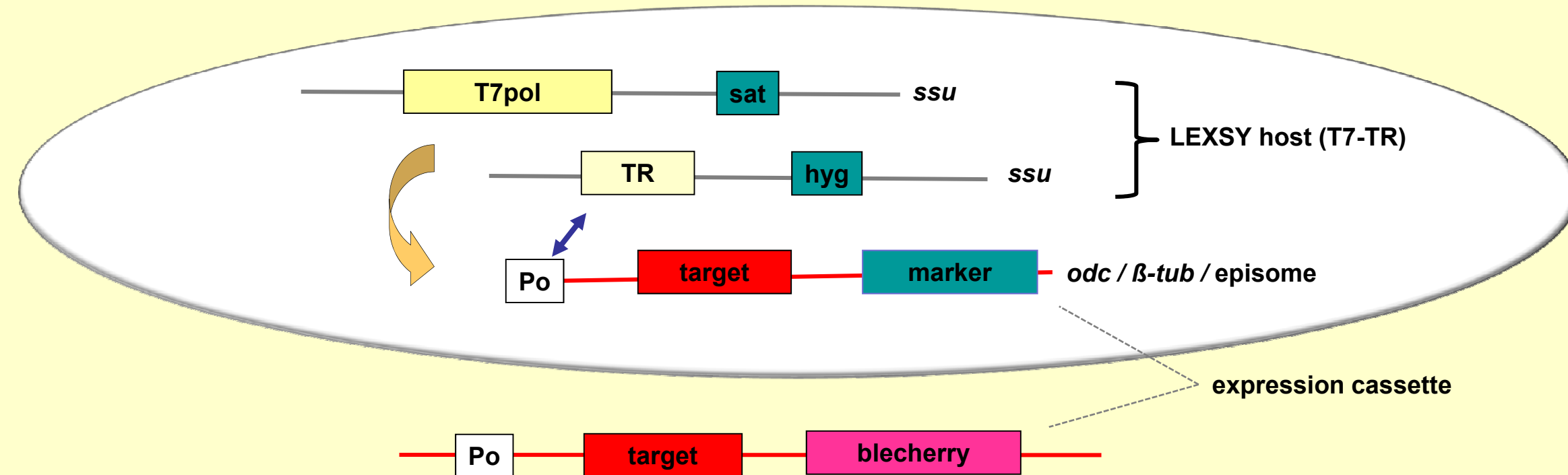
## Two configurations of *in vivo* LEXSY

### Constitutive expression of target genes



Transcription by strong host RNA polymerase I. Target genes are flanked by host-specific calmodulin (*cam*) splicing signals. Due to high homology of *ssu* genes used for integration, LEXSY vectors are also functional in other *Leishmania* sp. as *L. amazonensis*, *L. donovani*, *L. infantum*, *L. major*, *L. mexicana*, and also in the plant parasite *Phytomonas serpens*. Four selection markers (LEXSY NTC, Hyg, Bleo and Neo) permit co-expression of multi-subunit proteins.

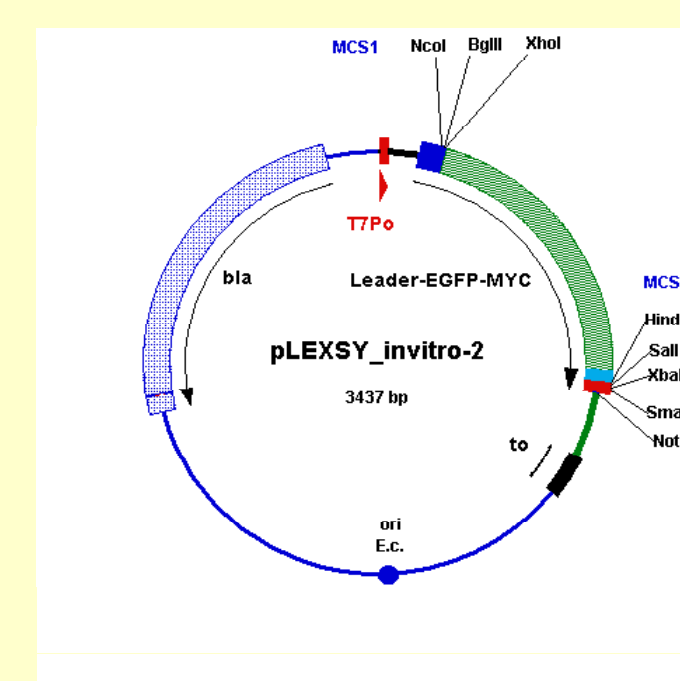
### Inducible expression of target genes



Uncoupling from LEXSY host regulation due to transcription by heterologous T7 RNA polymerase controlled by TET repressor. Selection with LEXSY Bleo or Neo. Blecherry marker permits selection and online monitoring of induction.

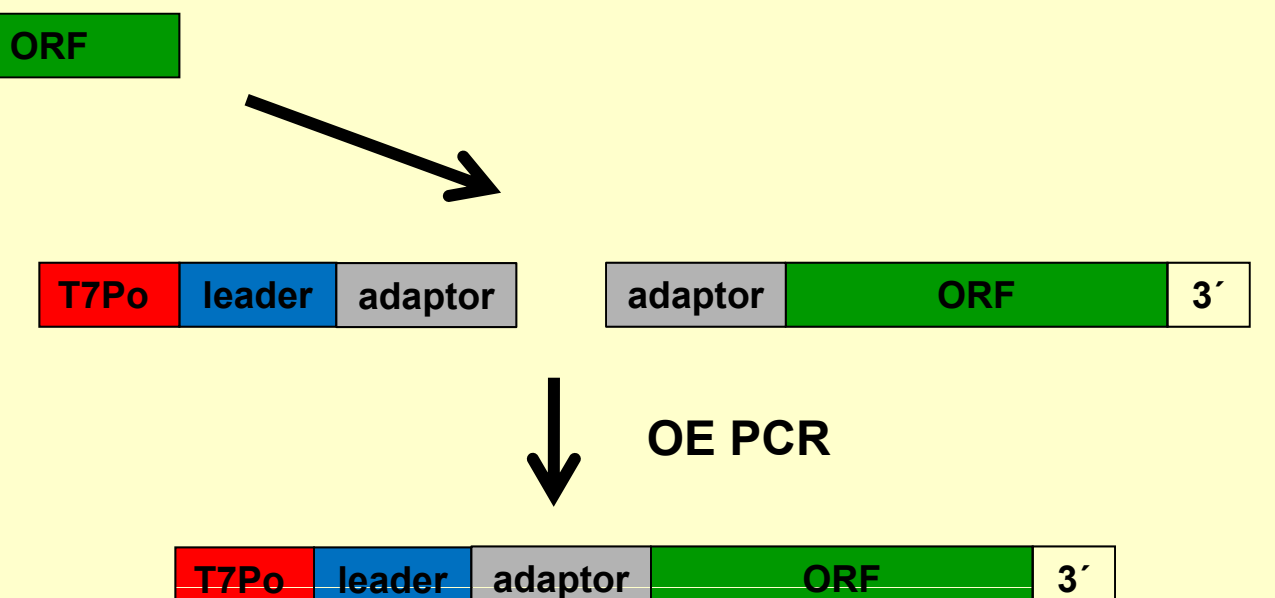
## Two configurations of *in vitro* LEXSY

### Plasmid based *in vitro* LEXSY - for high yields -



Template generation by cloning of target genes into plasmid vector

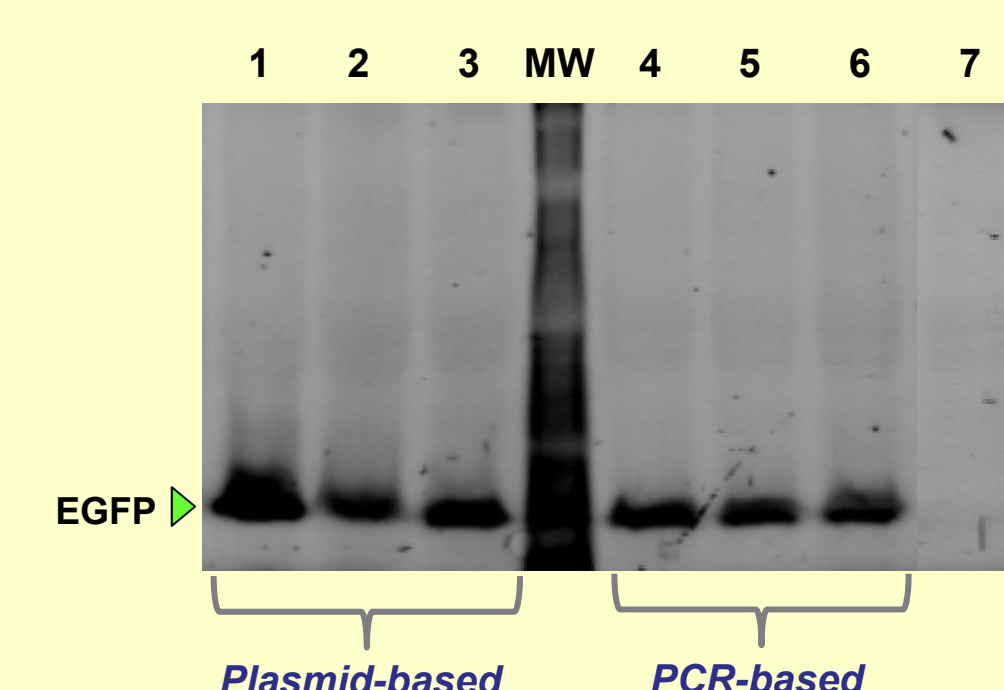
### PCR based *in vitro* LEXSY - for high throughput -



Template generation by direct PCR amplification of target DNA by overlap extension (OE) PCR



Cell-free production of proteins in transcription-translation coupled LEXSY cell extracts



Cell-free production of EGFP reference protein with plasmid-based (lanes 1-3) and PCR-based (lanes 4-6) *in vitro* LEXSY. Lane 7 negative control without template, MW molecular size marker. The *in vitro* reactions were carried out for 2 h at 20°C, resolved on 12% SDS-PAGE and EGFP was *in situ* visualized on a UV transilluminator.

## Selected applications of LEXSY in *Kinetoplastida* research



Further selected references:  
Breitling *et al.* (2002) Non-pathogenic trypanosomatid protozoa as a platform for protein research and production. *Protein Expression and Purification* 25: 209-218

Kushnir *et al.* (2005) Development of an inducible protein expression system based on the protozoan host *Leishmania tarentolae*. *Protein Expression and Purification* 42: 37-46

Lukeš J *et al.* (2006) Translational initiation in *Leishmania tarentolae* and *Phytomonas serpens* (Kinetoplastida) is strongly influenced by pre-ATG triplet and its 5' sequence context. *Molecular & Biochemical Parasitology* 148: 125-32

Mureev *et al.* (2009) Species-independent translational leaders facilitate cell-free expression. *Nature Biotechnology* 27: 747

Kovtun *et al.* (2010) Towards the Construction of Expressed Proteomes Using a *Leishmania tarentolae* Based Cell-Free Expression System. *PLOS one* 5: e14388

Kushnir *et al.* (2011) Artificial linear episome-based protein expression system for protozoan *Leishmania tarentolae*. *Molecular & Biochemical Parasitology* 176: 69

Jena Bioscience GmbH  
Loebstedter Str. 80  
07749 Jena, Germany  
Tel.: +49-3641-628-5000  
Fax: +49-3641-628-5100  
e-Mail: [info@jenabioscience.com](mailto:info@jenabioscience.com)  
[www.jenabioscience.com](http://www.jenabioscience.com)