

# FARM ANIMAL PROTEOMICS

KOŠICE, SLOVAKIA, 25-26 April 2013 *3rd Meeting*



## The LEXSY platform for recombinant protein expression

Reinhard Breitling

Jena Bioscience GmbH  
Loebstedter Str. 80  
07749 Jena, Germany  
Tel.: +49-3641-628-5000  
Fax: +49-3641-628-5100

e-Mail: [reinhard.breitling@jenabioscience.com](mailto:reinhard.breitling@jenabioscience.com)

# Shortcomings of conventional expression systems require alternative solutions



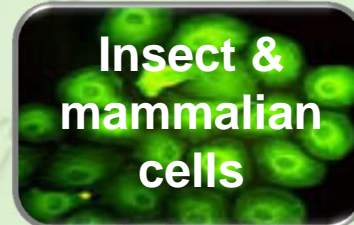
**Bacteria**

- Insufficient folding of complex proteins of higher organisms
- Inclusion bodies
- Lack of post-translational modifications



**Yeast**

- Posttranslational modifications differ largely from mammalian cells (high mannose)
- Problematic cell disruption



**Insect & mammalian cells**

- Laborious construction of over-expressing strains
- Expensive media
- Low growth rates
- Difficult scale-up



**Transgenic plants & animals**

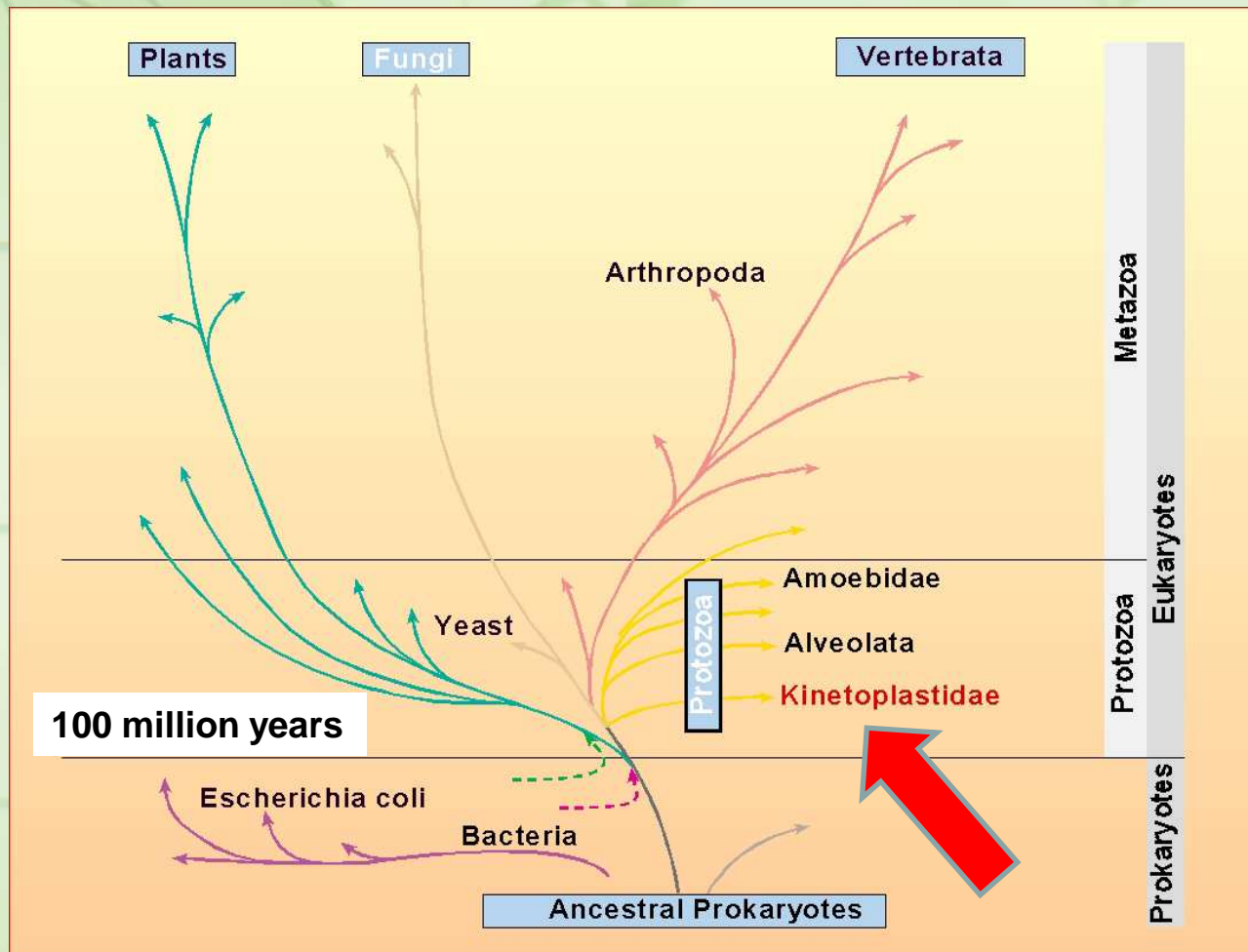
- Long developmental cycles
- Complex downstream processing
- Contamination problems (viruses)



**We need an expression system  
with eukaryotic machinery but bacterial robustness**



# Protozoa link the pro- and eukaryotic world



- Fully eukaryotic protein folding and modification
- molecular genetic manipulation developed
- Genome sequencing initiatives



Promising candidates among *Kinetoplastidae*

## *Leishmania tarentolae* was chosen as expression host

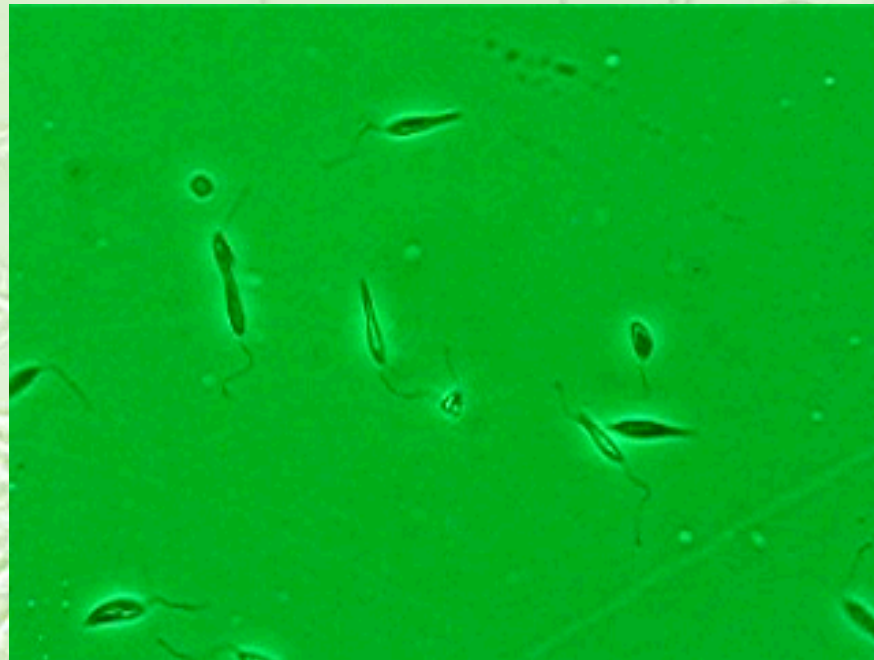


Host: *Tarentola mauritanica*

- unicellular flagellated protozoa
- parasite of lizard (*Sauroleishmania*)
- not pathogenic for mammals
- **biosafety group 1 organism**
- can be easily cultivated *in vitro*
- genome sequence published 2012

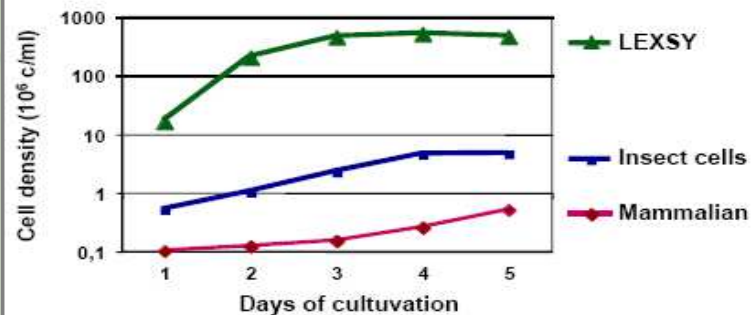
**“LEXSY”**

(Leishmania Expression  
System)

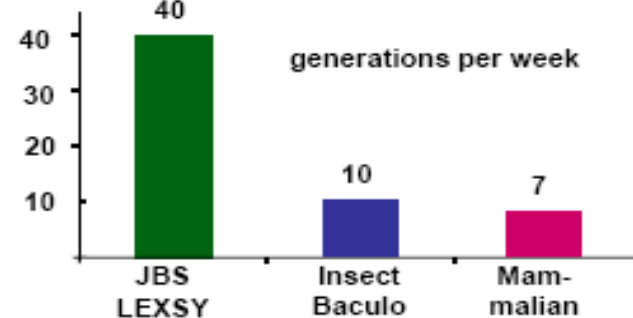


# LEXSY combines eukaryotic cellular machinery with bacterial robustness

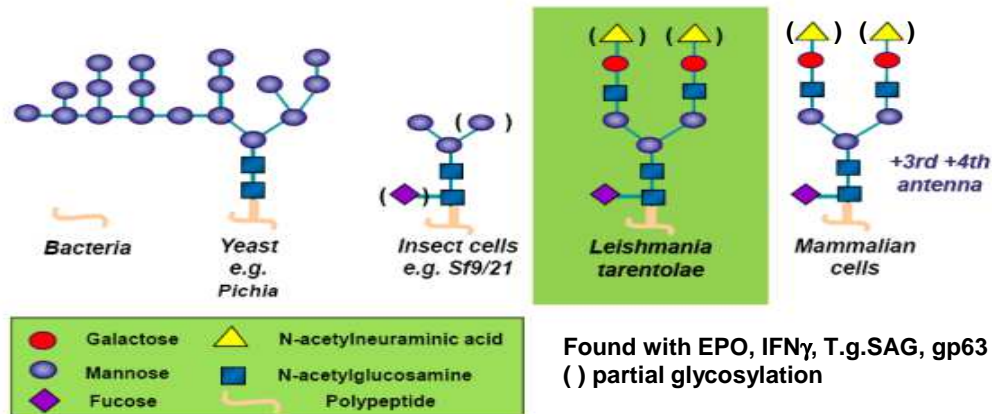
## LEXSY grows to higher densities than insect and mammalian cells



## LEXSY grows faster than insect and mammalian cells

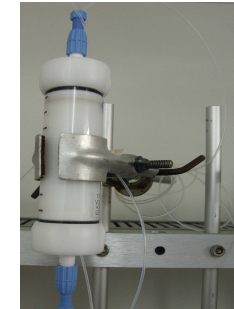
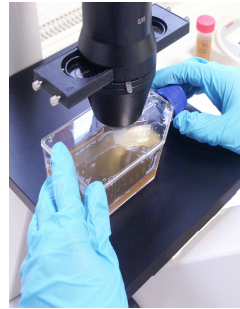
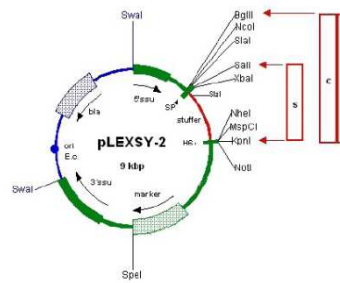


## LEXSY performs mammalian-type glycosylation





# With LEXSY - in six weeks from gene to protein



gene

**Expression  
plasmid  
construction**

**Transfection  
& selection**

**Expression  
evaluation**

**Scale-up**

**Protein  
purification**

protein

**Easy cloning  
in *E. coli***

**Versatile  
LEXSY  
expression  
vectors**

**1 week**

**Reliable  
electropora-  
tion  
protocols**

**Clonal or  
polyclonal  
selection**

**2 weeks**

**Constitutive  
or inducible  
  
Intracellular  
or secretory**

**Fluoresc.  
monitoring**

**1 week**

**Fully  
adapted to  
common  
fermentation  
technology**

**Up to 100  
litres tested**

**1 week**

**One-step  
affinity  
purification**

**Conventional  
techniques**

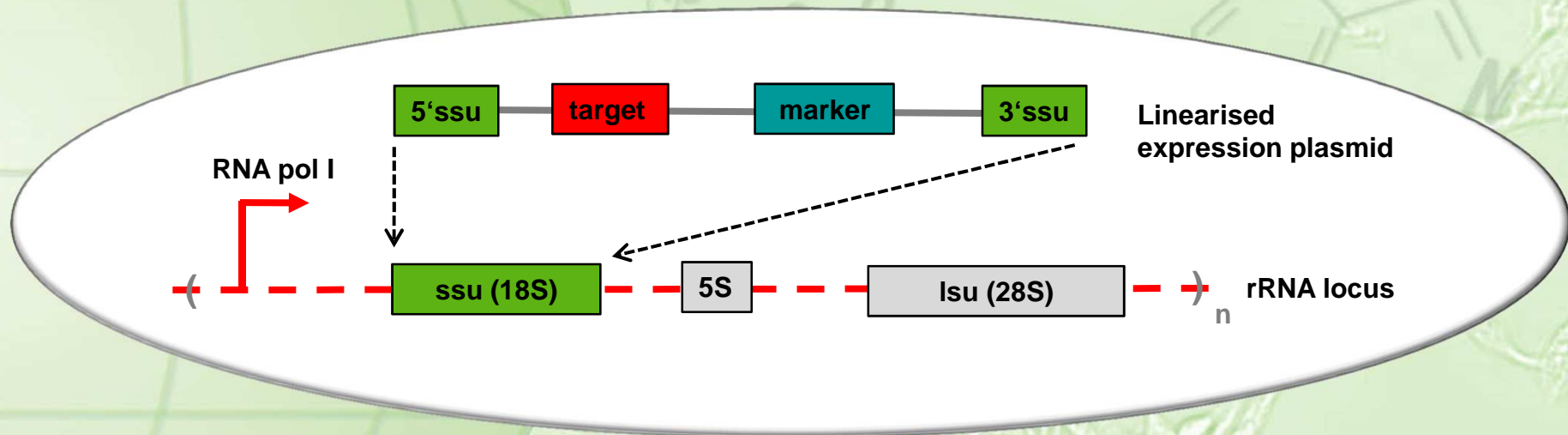
**1 week**



**LEXSY enables short evaluation cycles**

# Constitutive expression is the standard application for most proteins

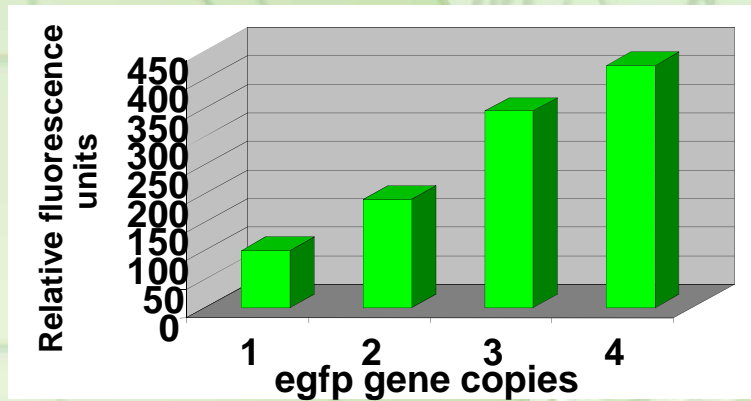
- *ssu* integration (multicopy 18S rRNA cluster)
- transcription by host RNA polymerase I



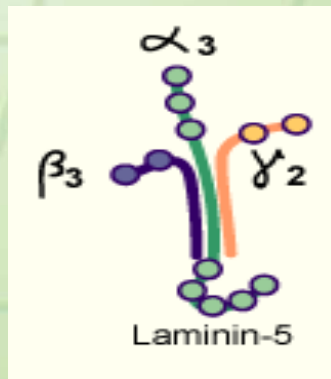
- RNA polymerase I: 10x stronger than RNA Polymerase II
- maximal activity during exponential growth (early product synthesis)
- four selection markers (LEXSY-NTC, -Bleo, -Hyg & -Neo)

# Co-integration of expression cassettes into *ssu* locus yielded additive effect of protein expression and allowed production of functional multi-subunit proteins

Expression levels



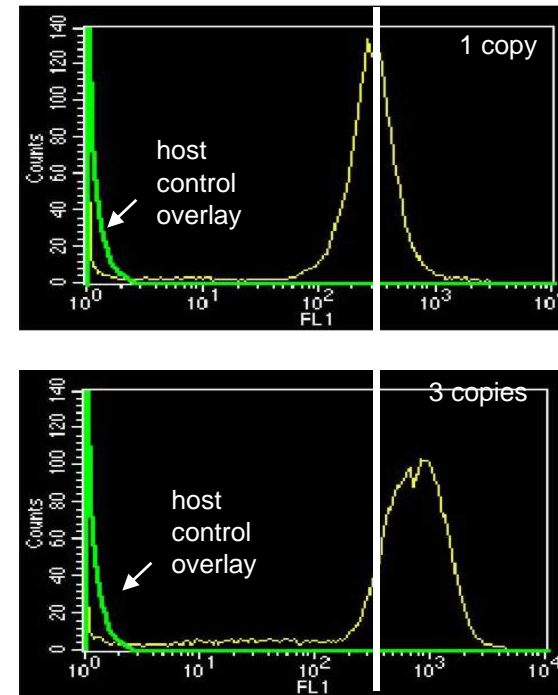
Also confirmed with eukaryotic target proteins EPO and Cu/Zn SOD



## Expression of multi-subunit proteins

Phan *et al.* (2009) **420 kDa** Laminin-332 heterotrimer

FACS analysis





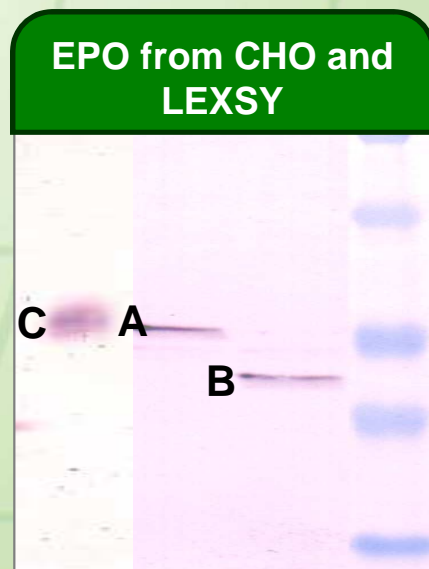
## Case study

# Human Erythropoietin was exceptionally homogeneously glycosylated in LEXSY

- Completely secreted to the culture medium
- Natively processed at the N-terminus
- Biologically fully active
- Exceptionally homogeneously & mammalian-type N-glycosylated  
(biantennary fully galactosylated  $\text{Man}_3\text{GlcNAc}_2\text{core-}\alpha\text{-1,6-fucosylated}$  structure)



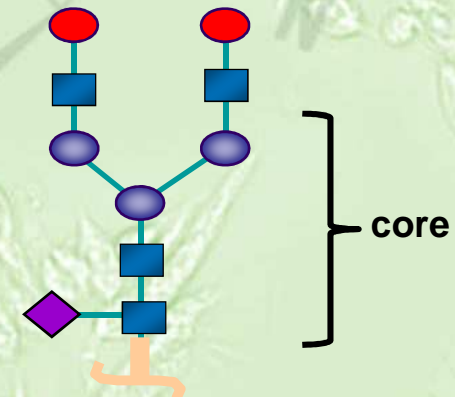
$1.2 \times 10^5$  U/mg



A: homogeneously glycosylated EPO from LEXSY

B: N-deglycosylated EPO from LEXSY

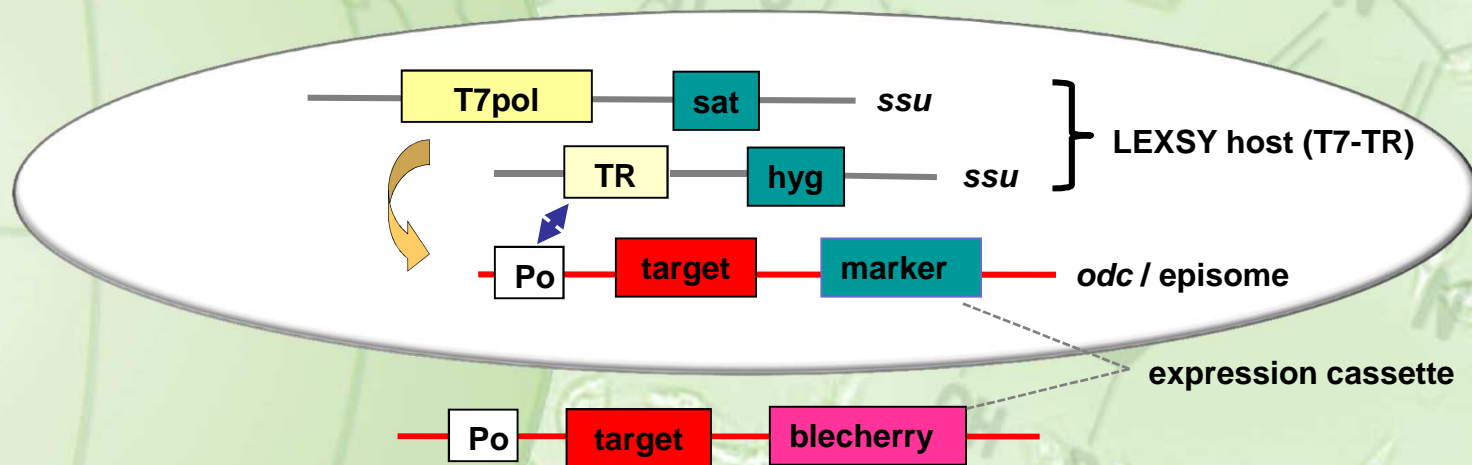
C: heterogeneously glycosylated EPO from CHO



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

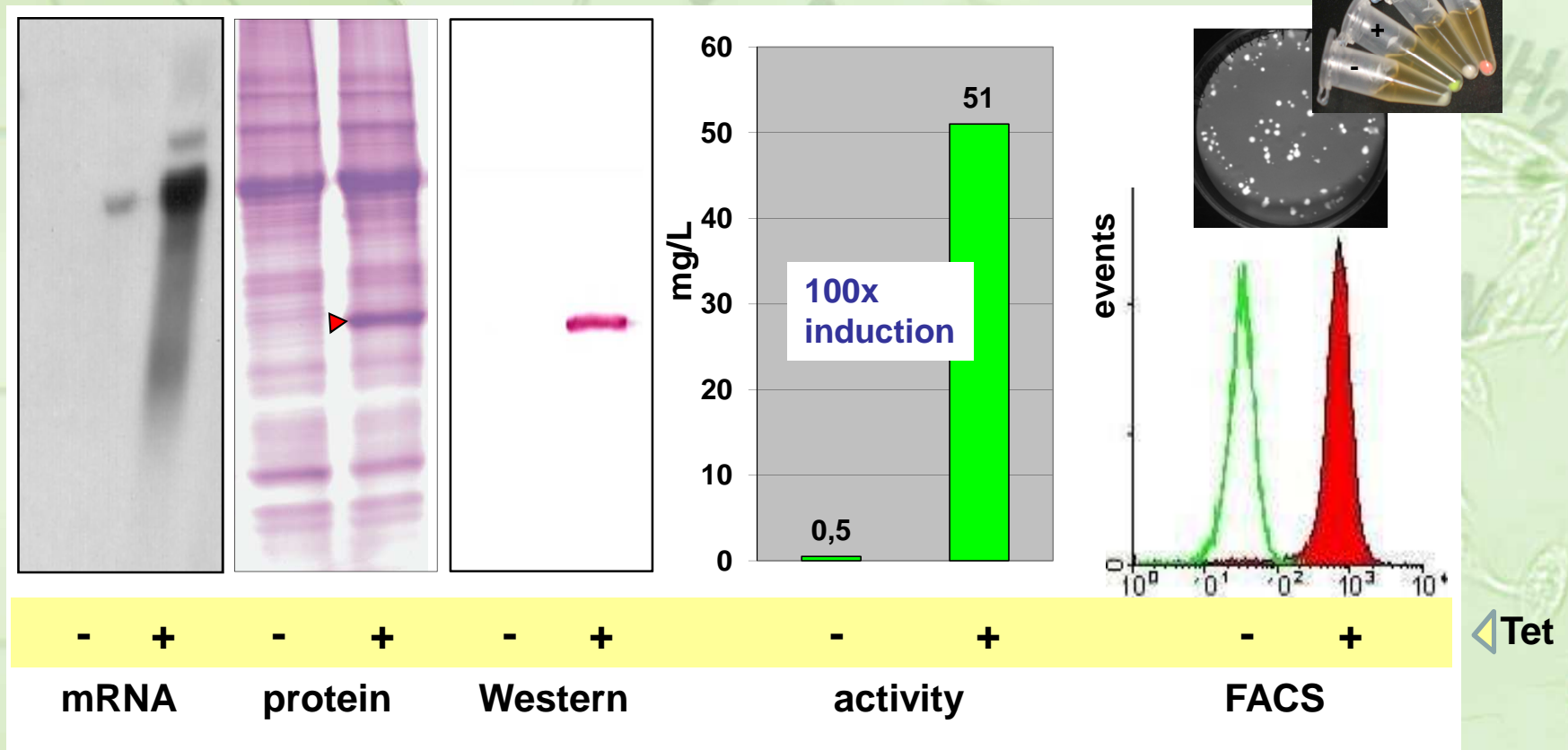
# Inducible expression is the application for difficult to express proteins

- *odc* integration or episomal propagation
- transcription by T7 RNA polymerase
- control by TET repressor



- uncoupling from cellular down-regulation
- much stronger than RNA Pol I
- tetracycline induction at any time during cultivation
- fluorescent marker (blecherry) for in-process monitoring

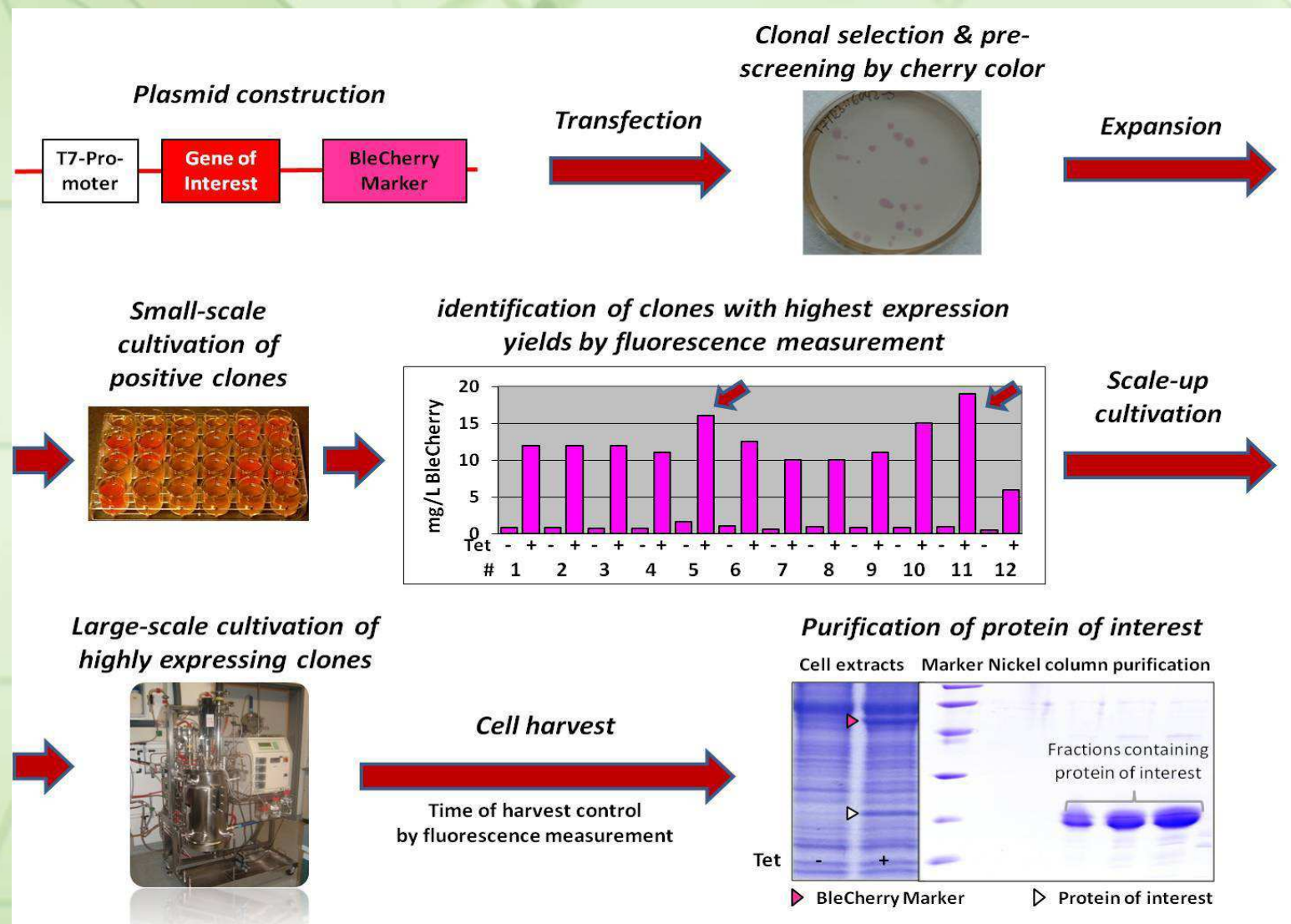
# Efficient expression control in inducible LEXSY from mRNA to protein



- induction profiles and yields stable over > 500 generations
- all cells induced



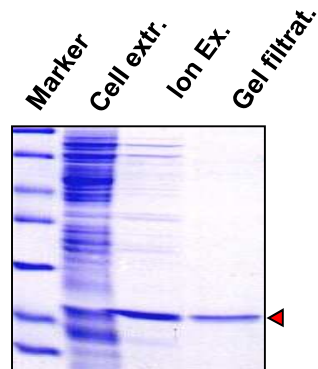
# LEXSY BleCherry architecture enables efficient screening and online monitoring of induction



# High-level protein expression in all four LEXSY formats

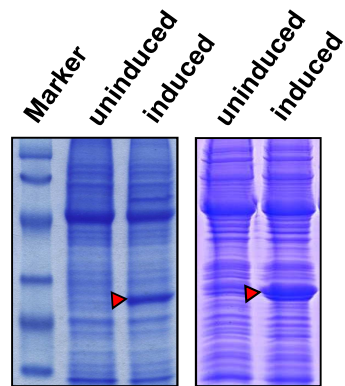
Constitutive

## Intracellular expression



Protein A

Two-step purification from lysed cells to purity > 95%

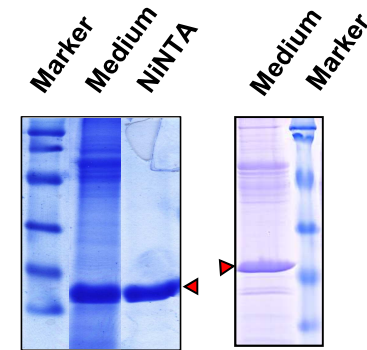


Protein D

Protein E

Up to 100 fold induction by addition of tetracycline

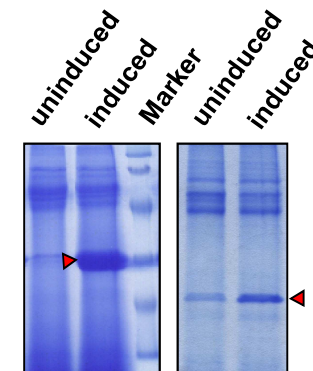
## Secretory expression



Protein B

Protein C

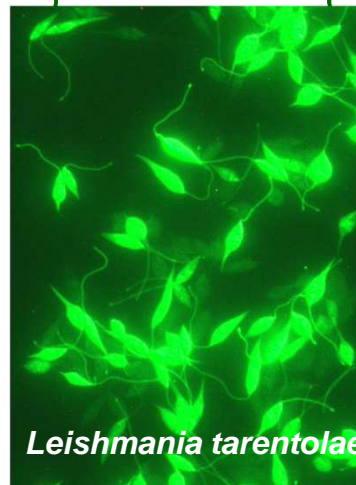
Protein of interest efficiently secreted to culture medium



Protein F

Protein G

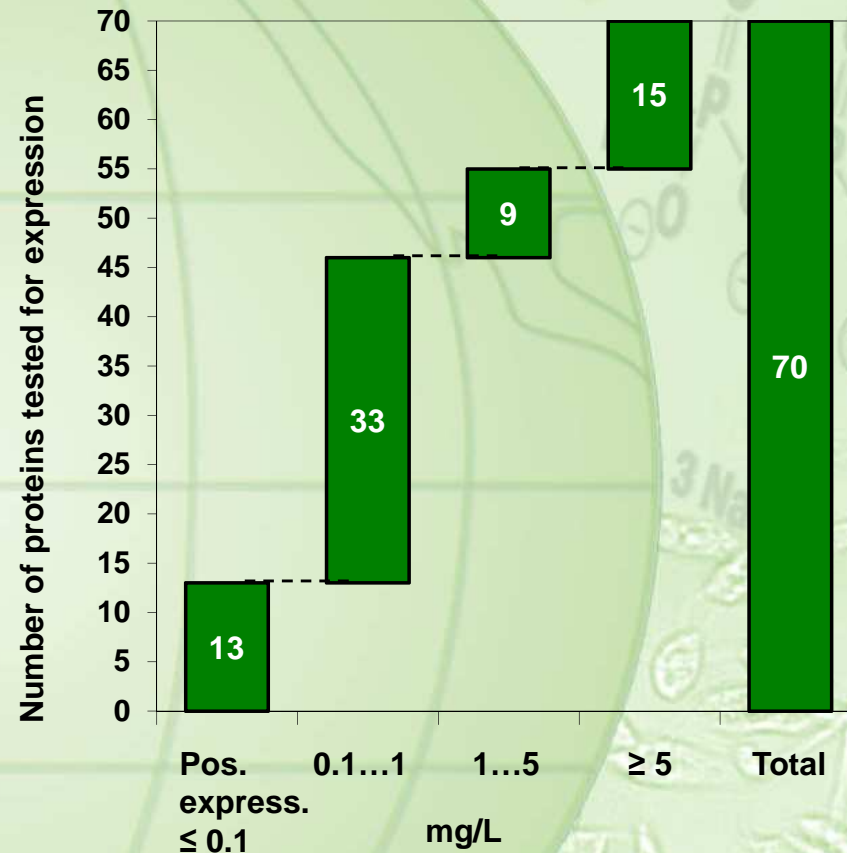
Upon Induction protein of interest becomes major component of cell culture medium



*Leishmania tarentolae*

► Protein of interest

**80% of target proteins expressed at > 0.1 mg/L**  
**1/3 > 1 mg/L, yields of up to 500 mg/L reached**



**40 cytosolic proteins**  
**24 secretory proteins**  
**6 membrane proteins**

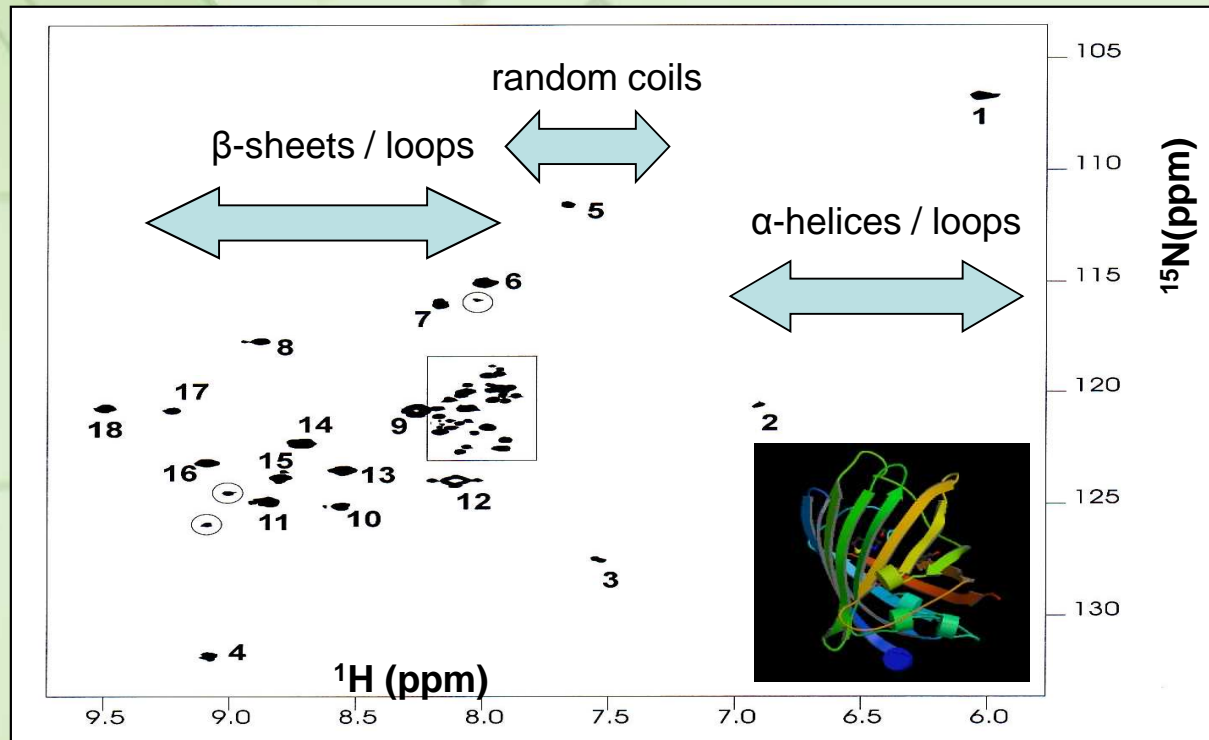
Target protein	Size kDa	Yield mg/L
Cytosplasmic proteins		
EGFP	28	300
SOD1	16	30
SPEE	35	30
p85 of PI3 kinase	85	3
smmyHC	154	1
Nuclear proteins		
T7 RNA Pol	100	1
Secreted proteins		
MHC II-β	30	500
CRP	23	44
SAG1&2	15/31	10
Fc fusion	39	10
MDP1	45	6
Laminin 332	420 (150+135+135)	0.5
Membrane proteins		
EGFP-Rab7 (mb-associated)	52	12
PDM9 (Type I)	43	0.5
BkrB2-GST (Type III TM7)	55	0.1



## Case study

## LEXSY enables *in vivo* protein labeling for NMR studies

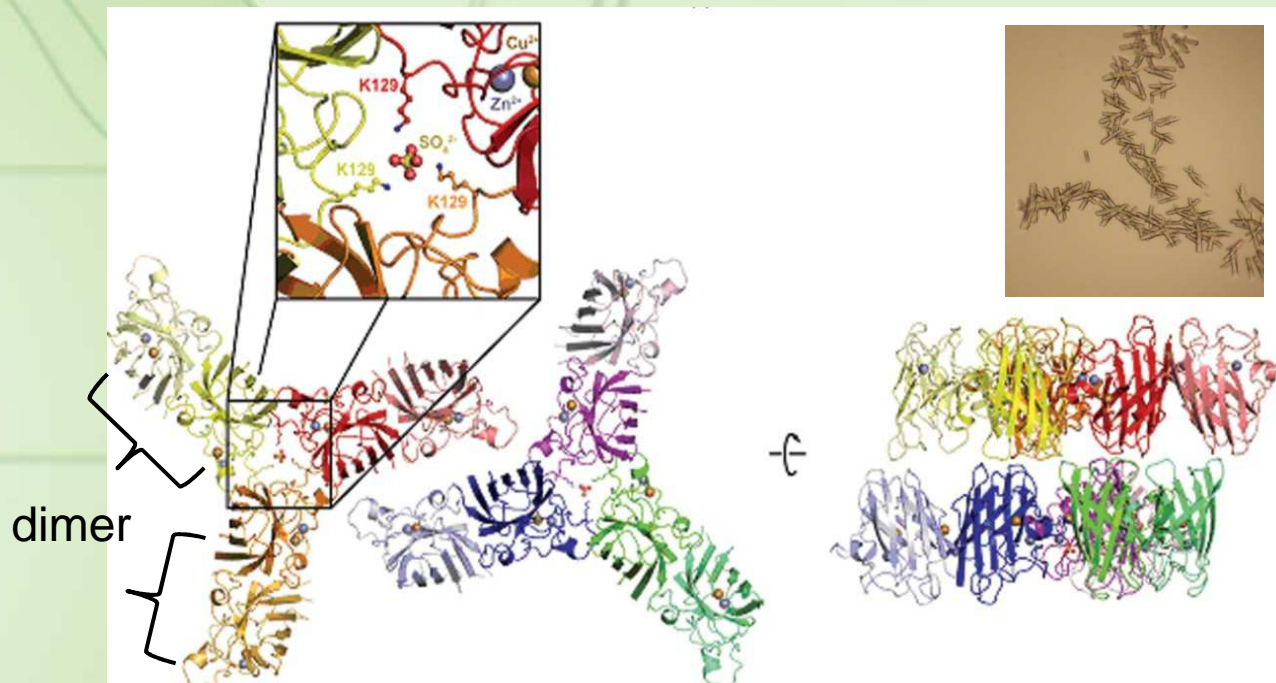
**- complete assignment of all 18  $^{15}\text{N}$ -Val residues in  $^{15}\text{N}$ -HSQC NMR of EGFP -**



- Niculae *et al.* (2006) Isotopic labeling of recombinant proteins expressed in the protozoan host *Leishmania tarentolae*. *Protein Expression and Purification* **48**: 167
- Foldynová-Trantírková *et al.* (2009) A Cost-effective Amino-acid-type Selective Isotope Labeling of Proteins Expressed in *Leishmania tarentolae*. *Journal of Biomolecular Structure & Dynamics* **26**: 755

**LEXSY is auxotrophic for 11 amino acids and can be grown in chemically defined media**

# Successful crystallography and X-ray analysis of LEXSY expressed proteins for structural biology



Structure determination of the **new P2<sub>1</sub>2<sub>1</sub>2<sub>1</sub> crystal form** of LEXSY-produced human Cu/Zn superoxide dismutase (SOD1). The asymmetric unit contains six SOD dimers arranged as two triangular wheels around sulfate ions. The wheels are arranged in a side-to-side fashion.

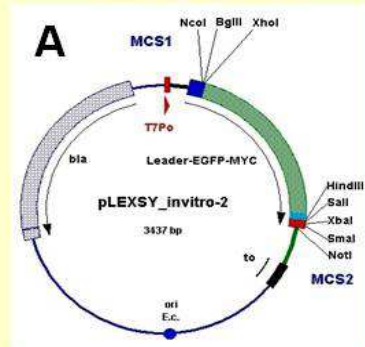
- Gazdag *et al.* (2010) Purification and crystallization of human Cu/Zn superoxide dismutase recombinantly produced in the protozoan *Leishmania tarentolae*. *Acta Crystallographica* **F66**: 871
- Dall & Brandstetter (2012) Activation of legumain involves proteolytic and conformational events, resulting in a context- and substrate-dependent activity profile. *Acta Crystallographica* **F68**: 24



# In Vitro LEXSY: Rapid cell-free protein production

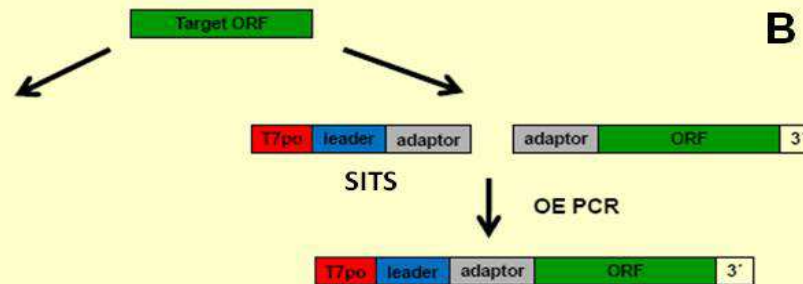
## Based on cell extracts of *L. tarentolae*

### 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 -

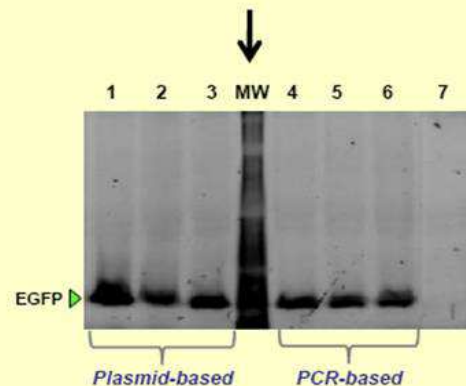
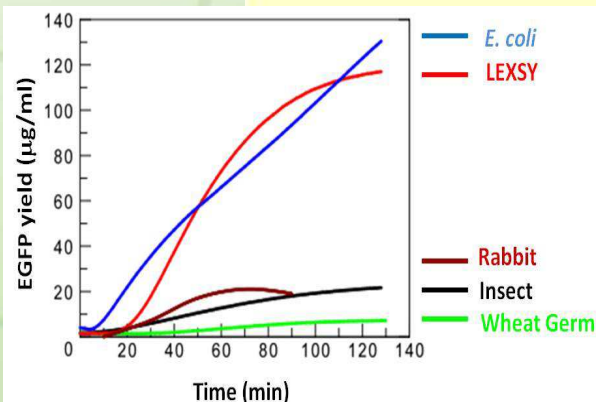


Template generation by direct PCR amplification of target DNA by overlap extension (OE) PCR

C



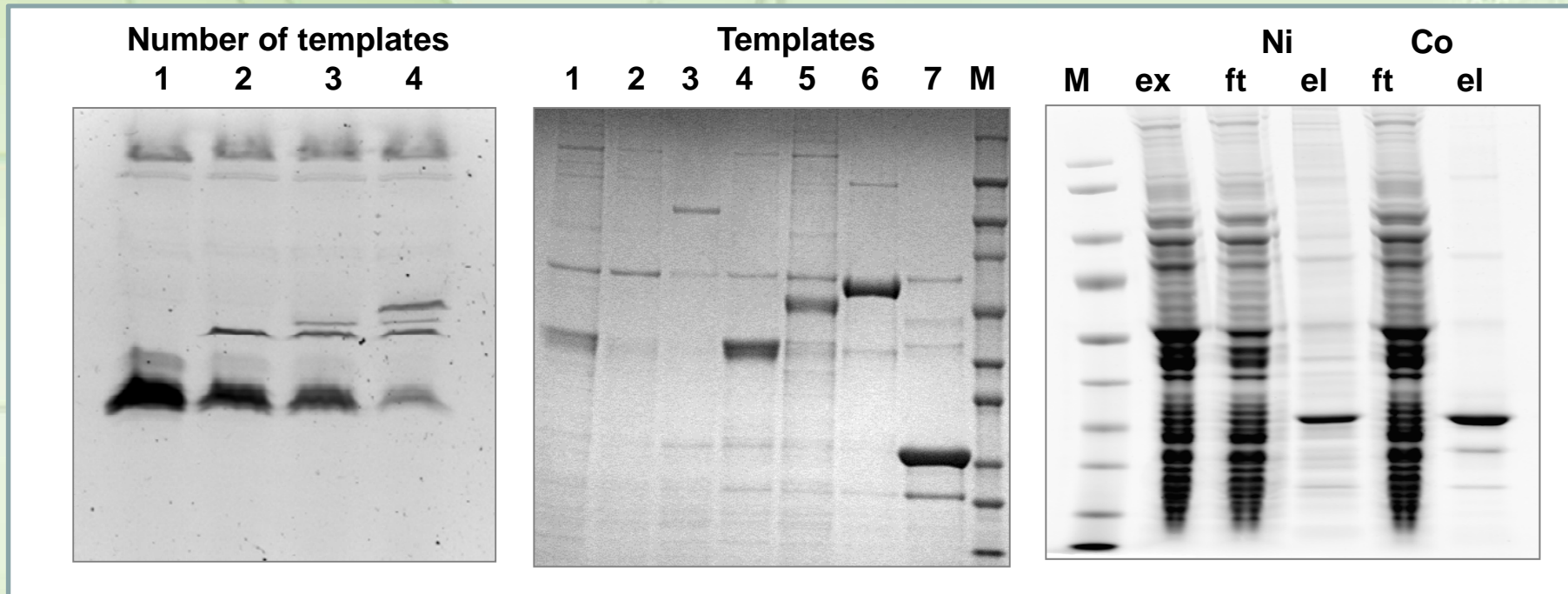
Cell-free production of proteins in transcription-translation coupled LEXSY cell extracts



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 *in situ* visualized on a UV transilluminator.



## ***In vitro* expressed EGFP fusion proteins can be visualised *in situ* and affinity purified**



In-gel ***in situ*** visualisation of EGFP fusion proteins **co-expressed** in the same extract

EGFP-**Cap matrix purification** of EGFP fusion proteins and detection by Coomassie staining

**Affinity purification** of *in vitro* synthesized EGFP by Co- and Ni-affinity matrices

- 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
- Kovtun *et al.* (2011) *Leishmania* cell-free protein expression System. *Methods* **55**: 58

## LEXSY - powerful protein expression system

- Robust fast growing unicellular host
- Eukaryotic protein synthesis / folding / modification
- Easy construction and rapid growth of LEXSY expression strains
- Short evaluation times
- Flexible expression solutions
- High yields and simple purification of target proteins
- LEXSY proteins for NMR and X-ray crystallography
- *In Vitro* LEXSY for rapid and parallel cell-free protein production

# Collaborators

## MPI molekul. Physiologie Dortmund

Kirill Alexandrov  
Susanna Kushnir  
Ion Cirstea  
Sergey Mureev  
Anca Niculae  
Mihai Gazdag  
Wulf Blankenfeldt  
Peter Bayer

## Univ. Queensland

Kirill Alexandrov  
Sergey Mureev  
Olexiy Kovtun

## FZMB Erfurt

Miriam Ebert  
Katrin Franke

## Washington Univ. St. Louis

Stephen M. Beverley

## University of Ghent

Nico Callewaert  
Roland Contreras

## Leibnitz Inst. Jena

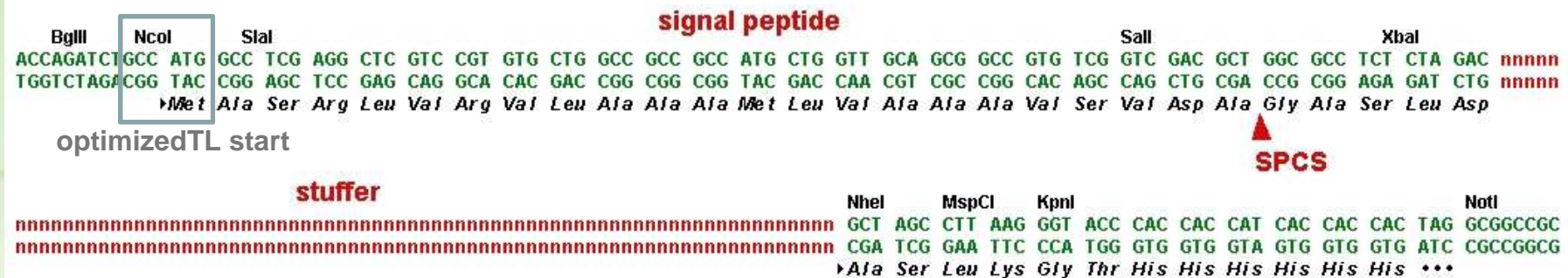
Karl-Heinz Gührs  
Alexander Tretjakov  
Vera Klujeva

## Univ. Applied Sciences Jena

Hans-Dieter Pohl  
Claudia Fritsche



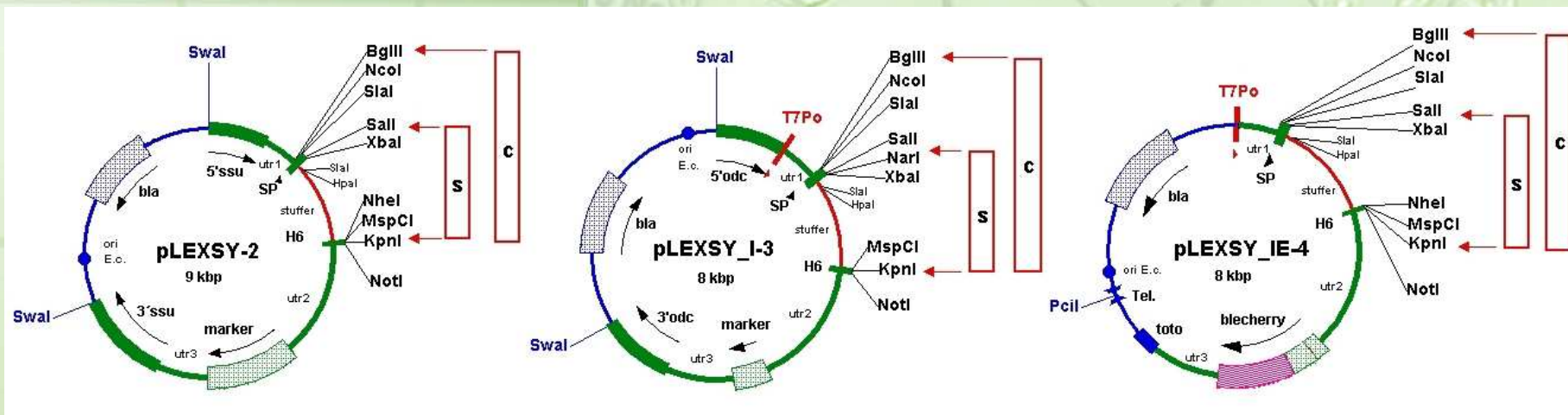
## Three LEXSY configurations for intracellular or secretory protein expression



## Constitutive integrative

## Inducible integrative

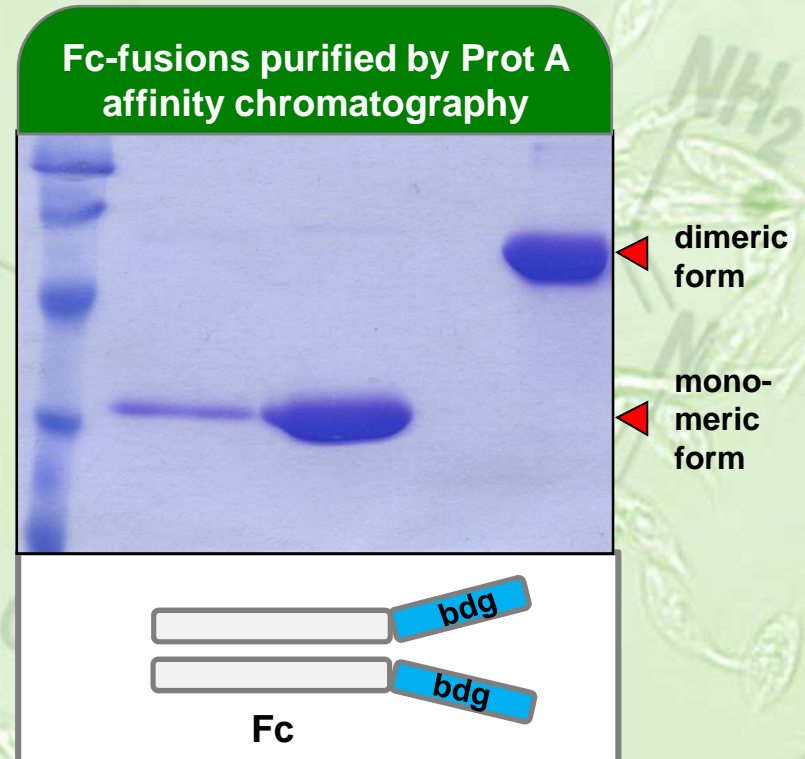
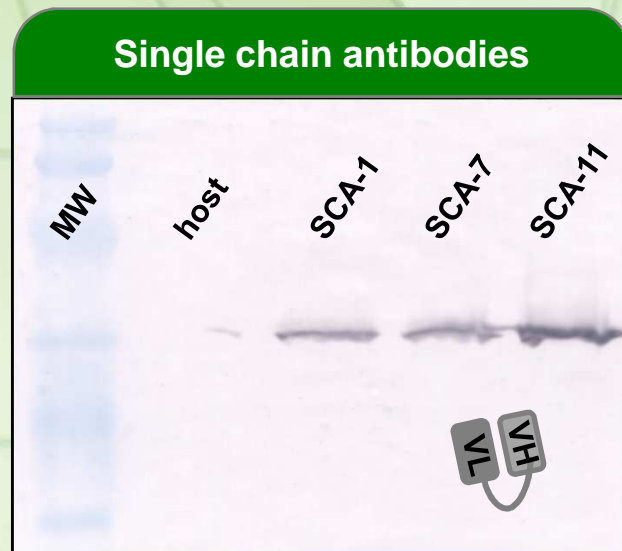
## Inducible episomal



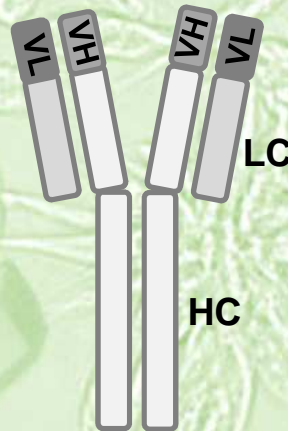
**C:** integration for cytosolic expression, **S:** integration for secretory expression  
**SP:** signal peptide of secreted acid phosphatase of *Leishmania mexicana*  
**SPC:** signal peptide cleavage site

## Case studies

# Different antibody constructs were expressed in LEXSY



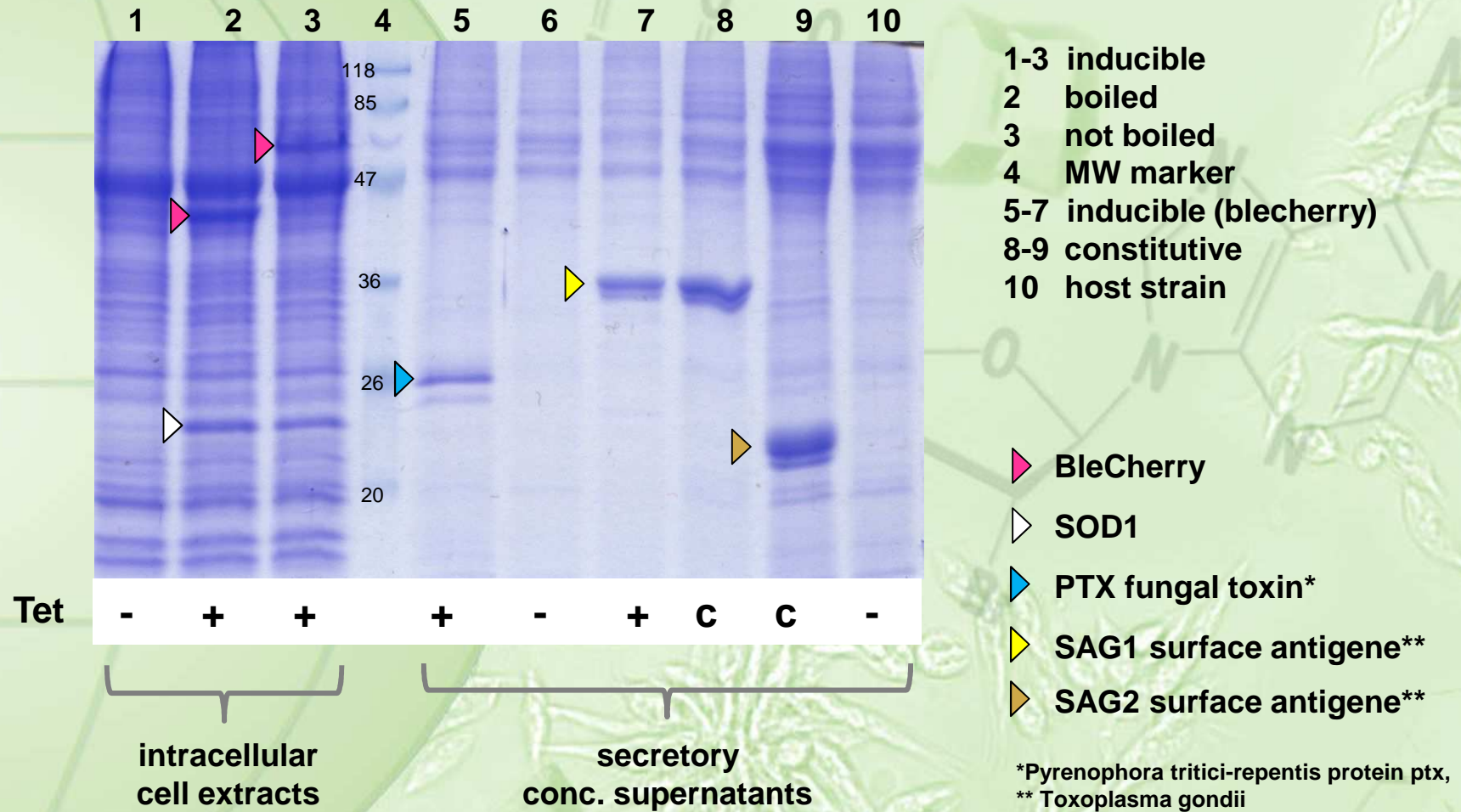
HC and LC of human IgG were expressed separately in different LEXSY strains





## Case studies

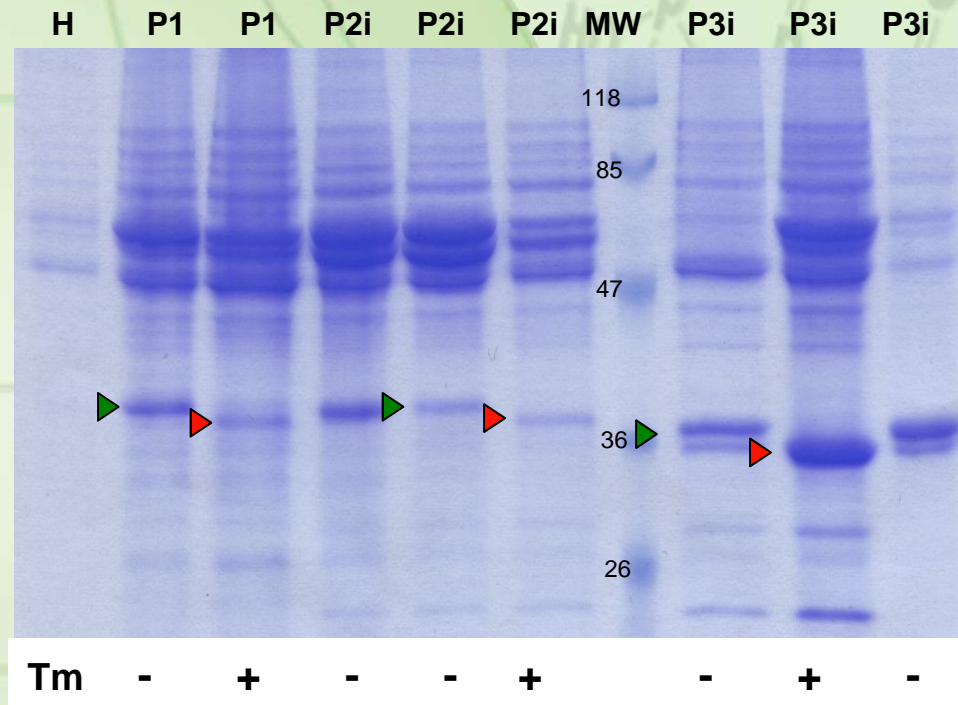
# Various target proteins were efficiently expressed in LEXSY





## Case studies

# Glycosylation of secretory target proteins was inhibited *in vivo* by addition of Tunicamycin



H LEXSY host (negative control)  
 MW prestained molecular size marker  
 P1 constitutive secretory expression  
 P2i inducible secretory expression  
 P3i inducible secretory expression (enzymatic deglycosylation was shown)  
 Tm Tunicamycin added to culture at 10 µg/ml

➤ Glycosylated target protein from LEXSY  
 ➤ Non-glycosylated target protein from LEXSY

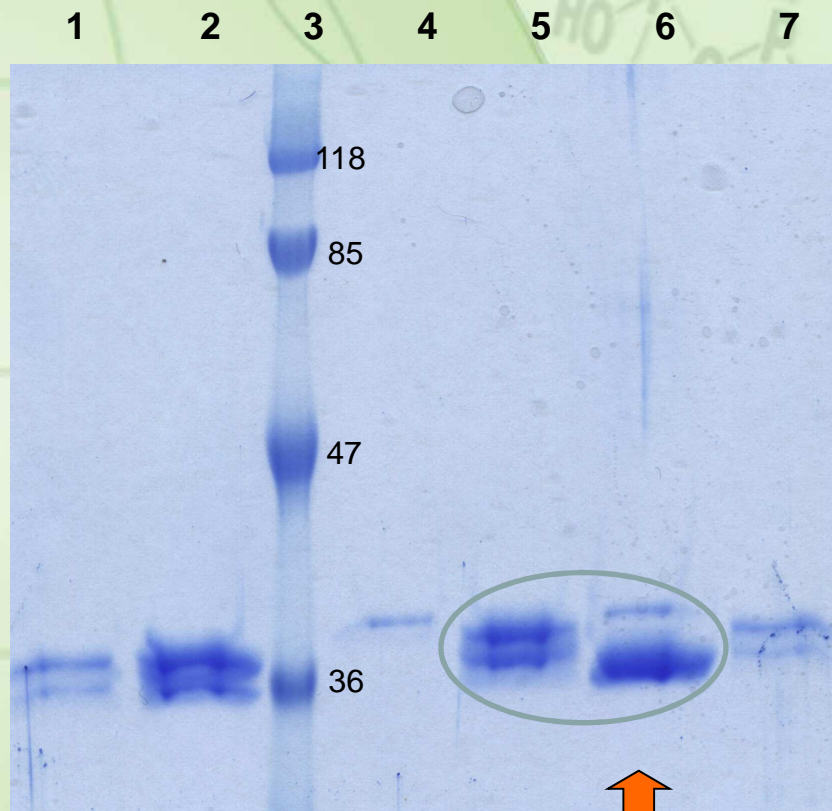
Concentrated culture supernatants of LEXSY clones

Mobility shift caused by addition of Tunicamycin to culture indicates glycosylation of target proteins in LEXSY

## Case studies

### LEXSY expressed protein was deglycosylated *in vitro*

Target protein was affinity-purified from culture supernatant and enzymatically deglycosylated



- 1 imidazol eluted POI
- 2 concentrated POI in PBS
- 3 prestained molecular size marker
- 4 N-Glycosidase F (138 ng = 250U) 36 kDa
- 5 POI in deglycosylation buffer
- 6 POI + N-Glycosidase F**
- 7 imidazol eluted POI

POI: target protein of interest

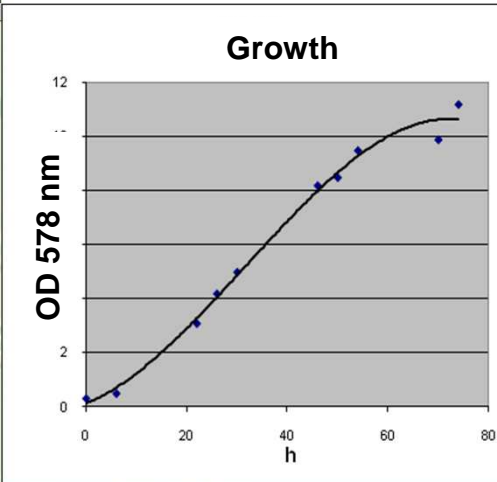
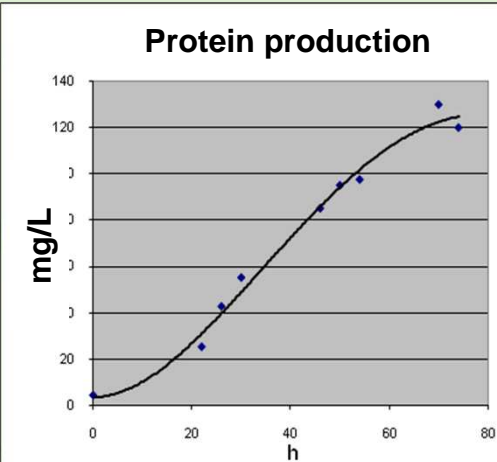
←  
← 2 glycoforms  
consistent with previous findings (EPO)

↑  
shift to one band



## Case studies

# LEXSY is compatible with common fermentation technology

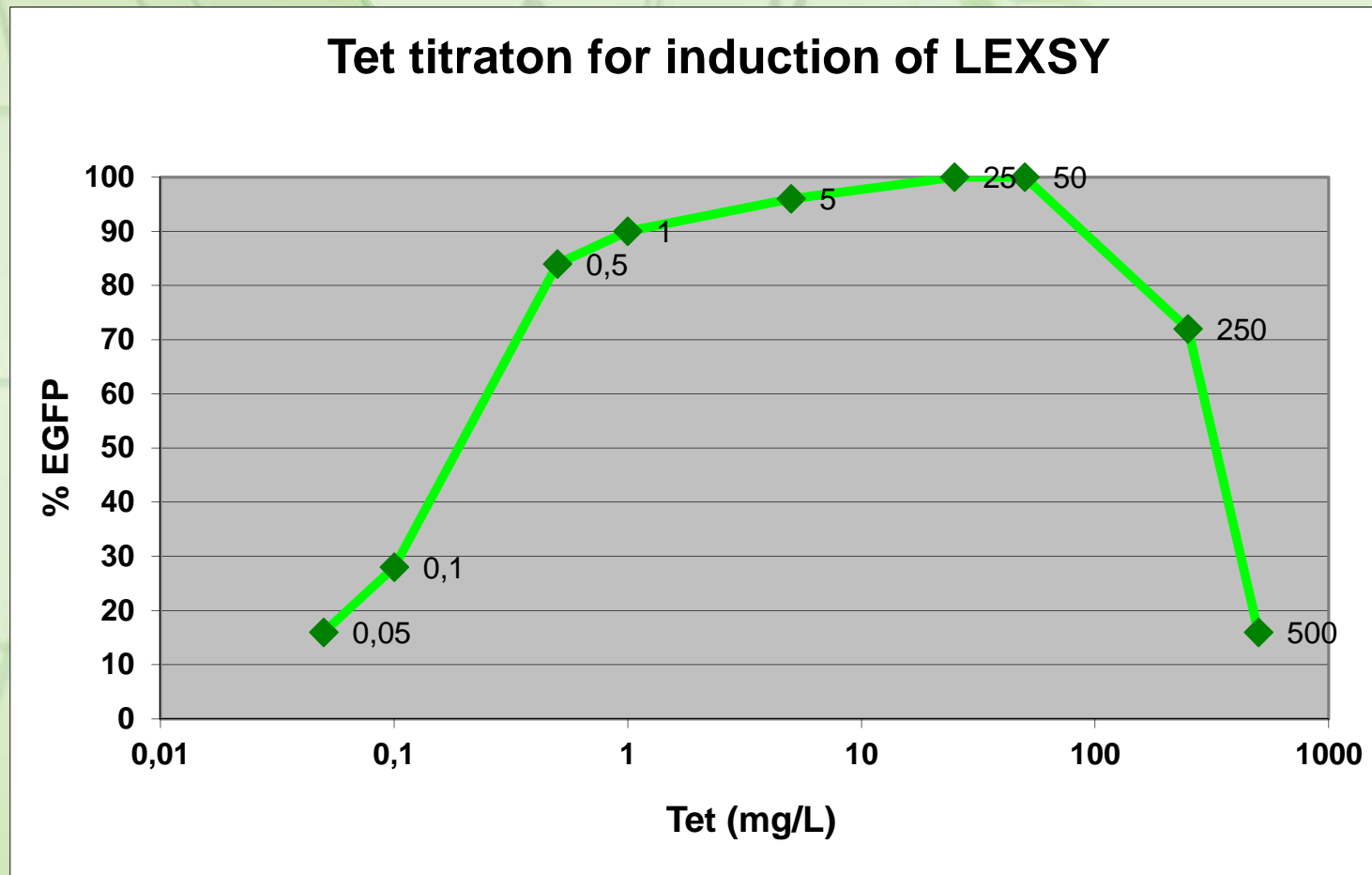


**Constitutive protein production parallels growth of expression culture  
>120 mg/L of culture reached at  $8 \times 10^8$  cells/ml**



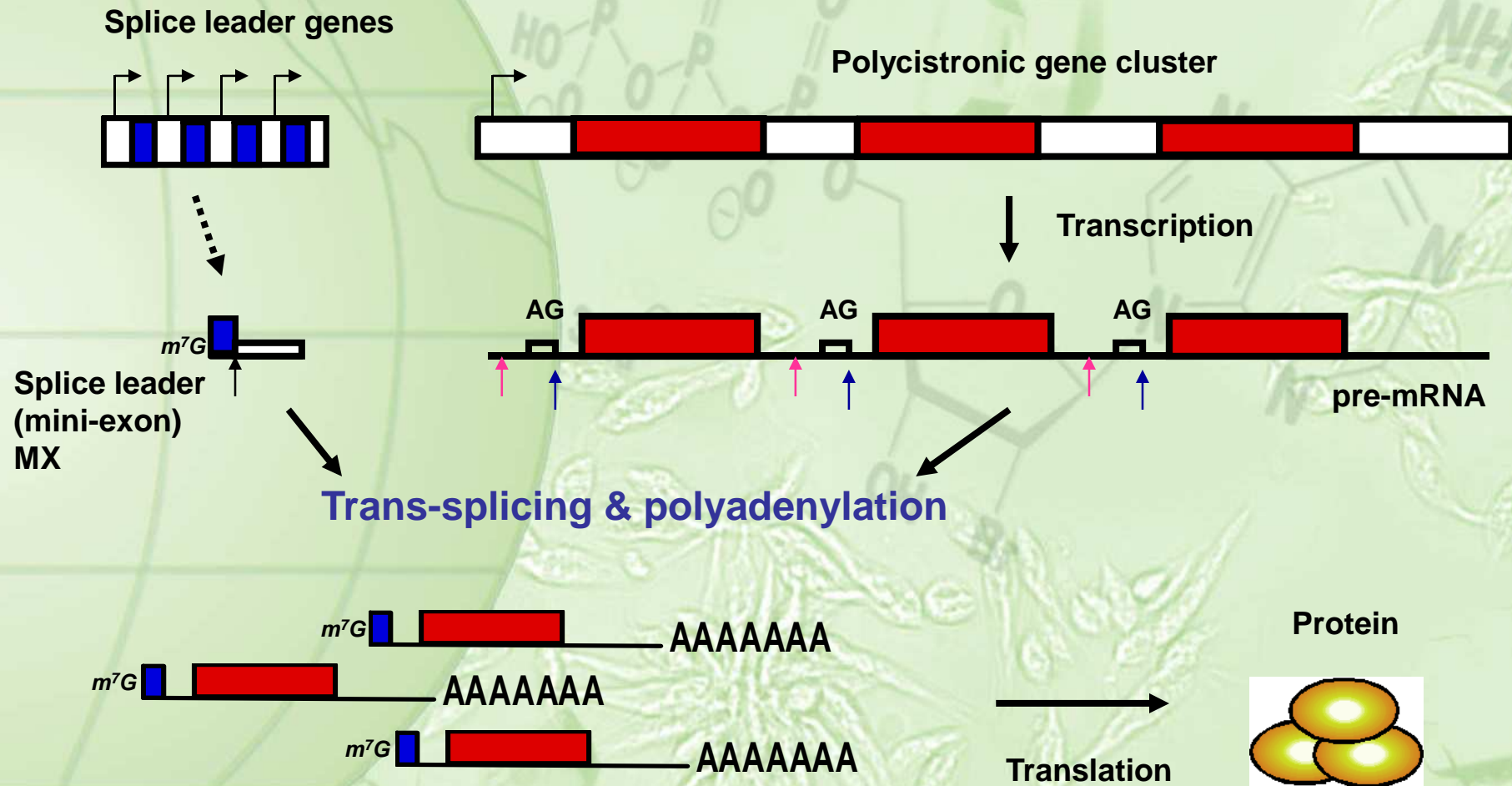
## Appendix

### Broad induction plateau in inducible LEXSY



## Appendix

# Transcription is uncoupled from RNA-processing in *Leishmania*

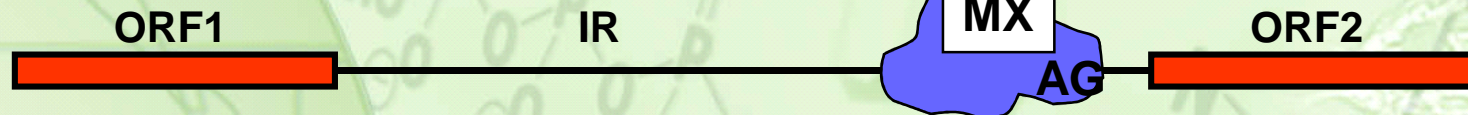




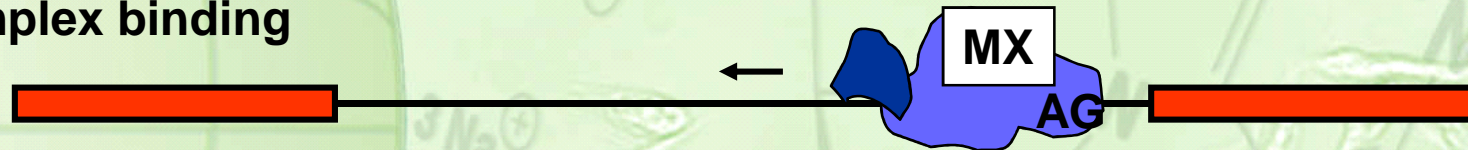
## Appendix

### Scanning model for coupling of splice site and poly(A) site selection in *Leishmania*

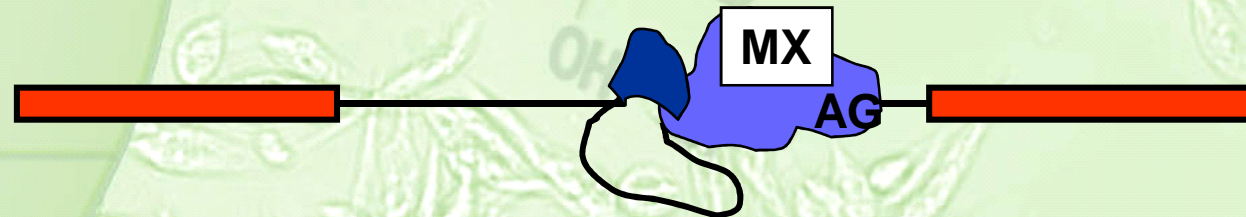
Spliceosome assembly



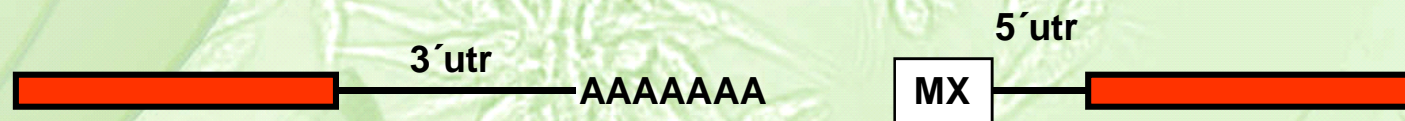
PolyA complex binding



Scanning 3' → 5'



Cleavages



MX: miniexon

After LeBowitz et al., 1993