



5th HALLE CONFERENCE

ON RECOMBINANT PROTEINS

19th – 20th
FEB 2015
LEOPOLDINA

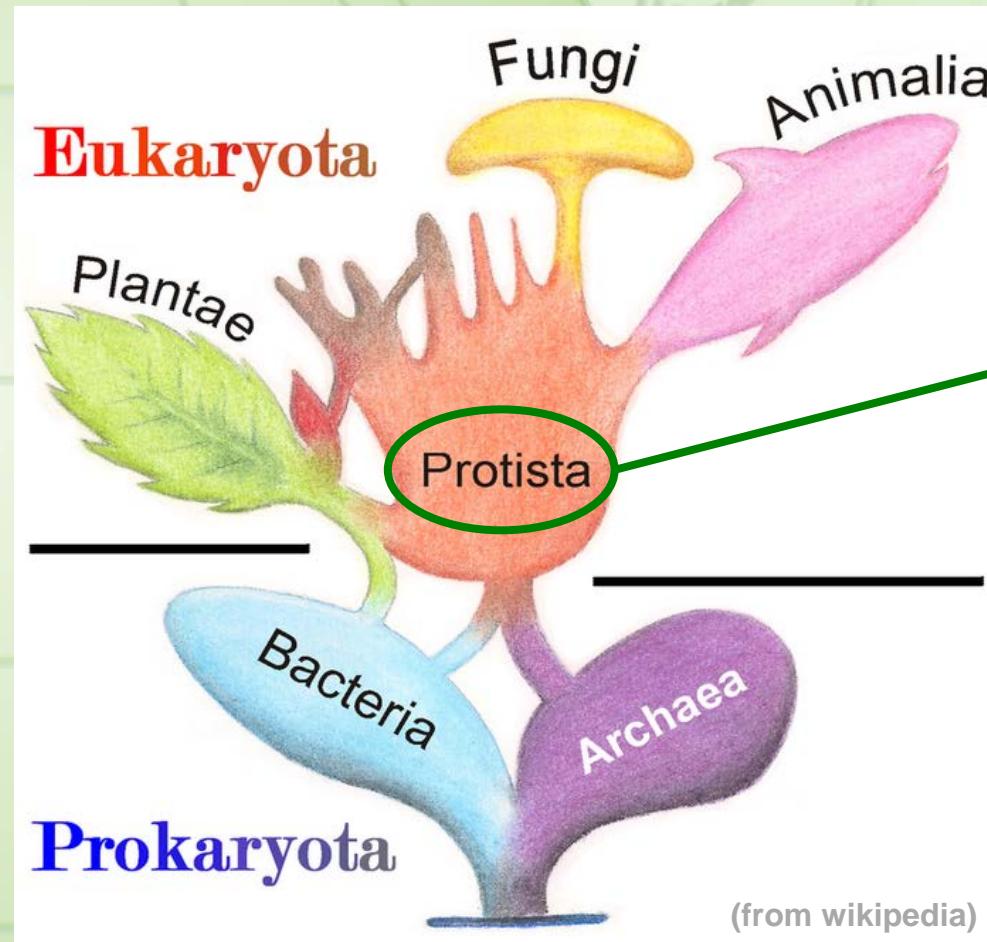
 Jena Bioscience

LEXSY taking off: Selected examples for protein production with *Leishmania tarentolae*

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

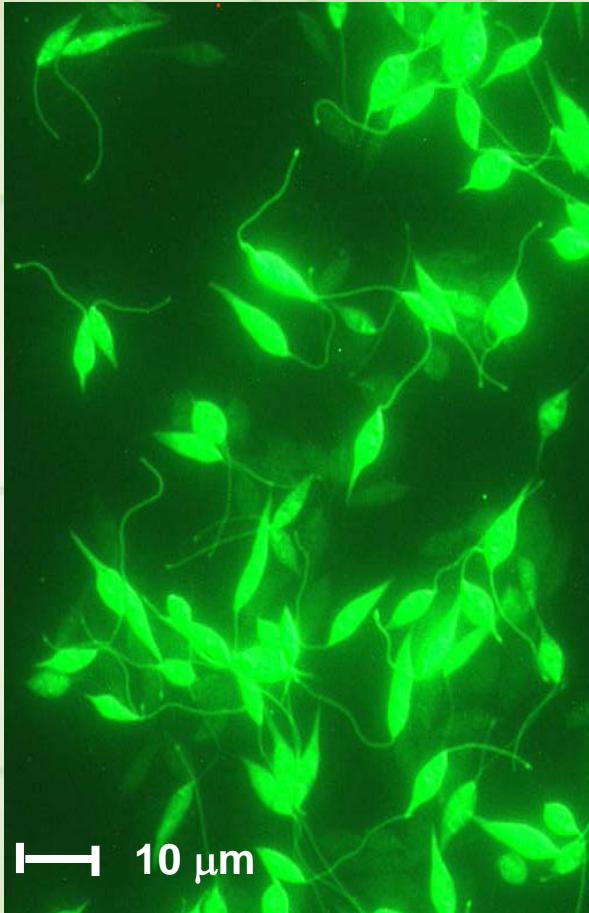
e-Mail: expression@jenabioscience.com
<http://www.jenabioscience.com/LEXSY>

Why not an expression system that **combines** prokaryotic easiness with eukaryotic capabilities?



- Large, diverse kingdom of organisms
- Many didn't look very promising...
- ...but one did:
Leishmania tarentolae

So *Leishmania tarentolae* became LEXSY



Leishmania tarentolae

- Nonpathogenic to mammals (S1-clearance)
- Very happy in culture (flasks and fermenters)
- Complete eukaryotic protein synthesis / folding / modification machinery (PTMs⁽¹⁾)

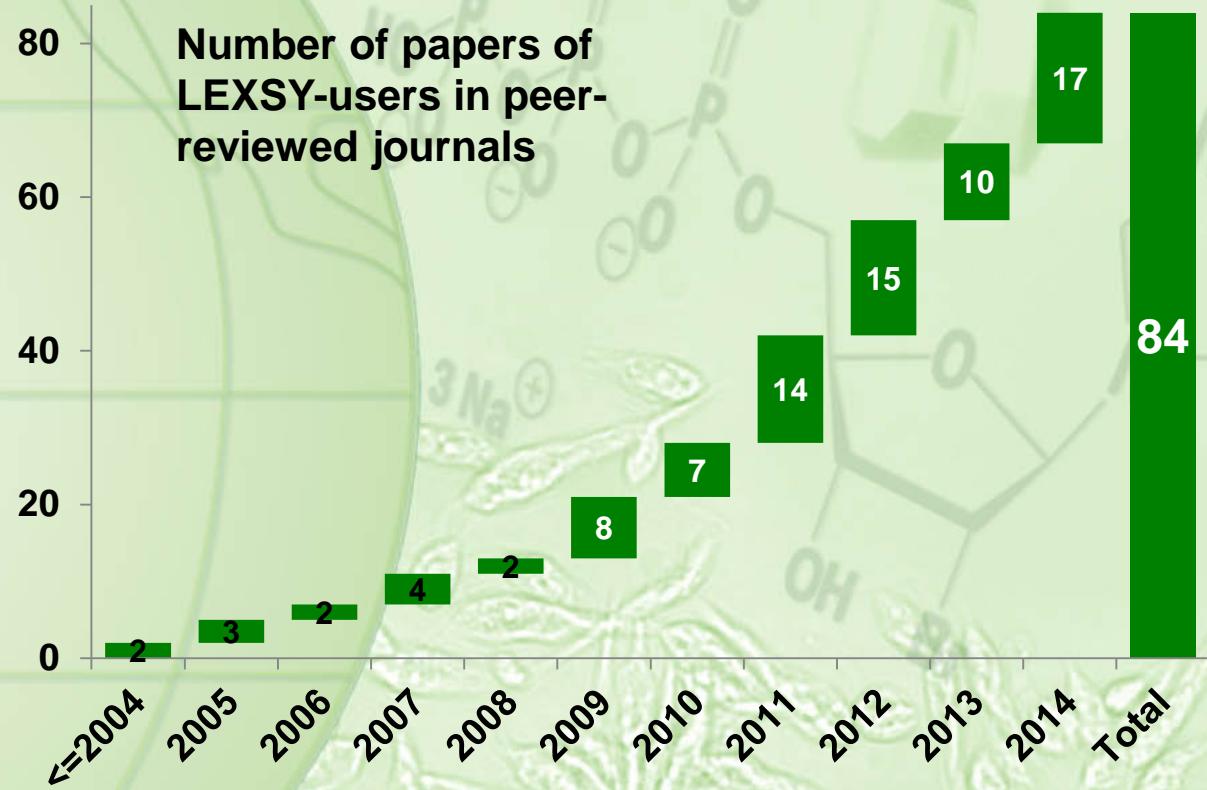
Not even 10 years of hard
lab work...

LEXSY

- Suitable for many *types* of proteins including membrane, cytosolic, nuclear and secreted proteins
- Approx. 80% positive expression projects with yields up to 500 mg per litre of culture
- Cell-free version (*in vitro* translation with LEXSY extracts)

(1) PTMs = post-translational modifications such as mammalian-type glycosylation, phosphorylation, glypiation (GPI anchoring), acetylation, prenylation, myristoylation, ADP-ribosylation, proteolytic processing, oligomerisation...

Soon more and more people started using LEXSY
To date, LEXSY is established in roughly 200 labs worldwide



Exponential take-off...? Last few minutes some selected LEXSY examples

LEXSY makes antigens, EPO, interleukin and interferon

Vaccines

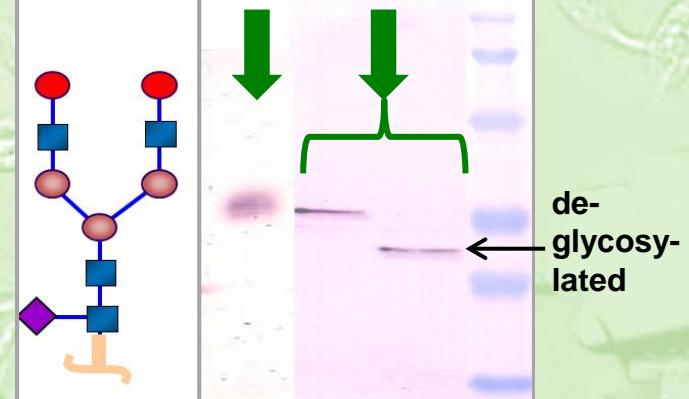
- Production of recombinant viral antigens (influenza⁽¹⁾, Hepatitis⁽²⁾, Papilloma Virus⁽³⁾)
- Production of recombinant parasite antigens (Leishmaniasis⁽⁴⁾)
- Whole LEXSY cells as live vaccine delivery vector in tumor models^(5,6)

Cytokines

- Human interferon gamma (IFN γ) with yields of 10 mg/L⁽⁸⁾
- N-glycosylated, antivirally active human interleukin 29 (IL-29⁽⁹⁾)
- Mammalian-type N-glycosylated homogeneous biologically active human EPO⁽⁷⁾

- (1) Pion et al. (2014) *Vaccine* 32: 5570
(2) Baechlein et al. (2013) *Journal of Virological Methods* 193: 238
(3) Hosseinzadeh et al. (2013) *Drug Delivery* 20: 190
(4) Chamakh-Ayari et al. (2014) *PLOS ONE* 9: e92708
(5) Saljoughian et al. (2013) *PLoS Neglected Tropical Dis.* 7: e2174
(6) Salehi et al. (2012) *Immunotherapy* 4: 1107
(7) Breitling et al. (2002) *Prot Expr and Purification* 25: 209
(8) Davoudi et al. (2011) *World J. Microbiol. Biotechnology* 27: 1893
(9) Taromchi et al. (2013) *Iranian Journal of Biotechnology* 11: 168

EPO from: CHO, LEXSY

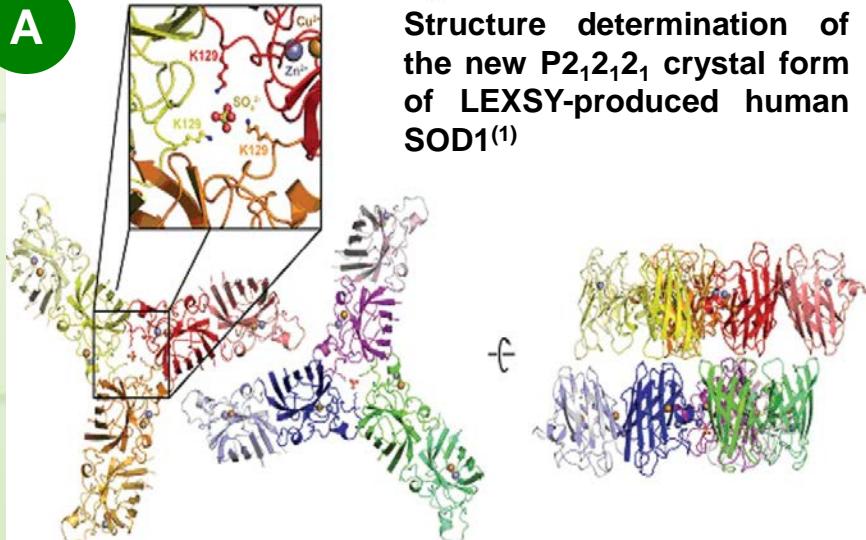


LEXSY-proteins can be crystallized and NMRed

Structural biology

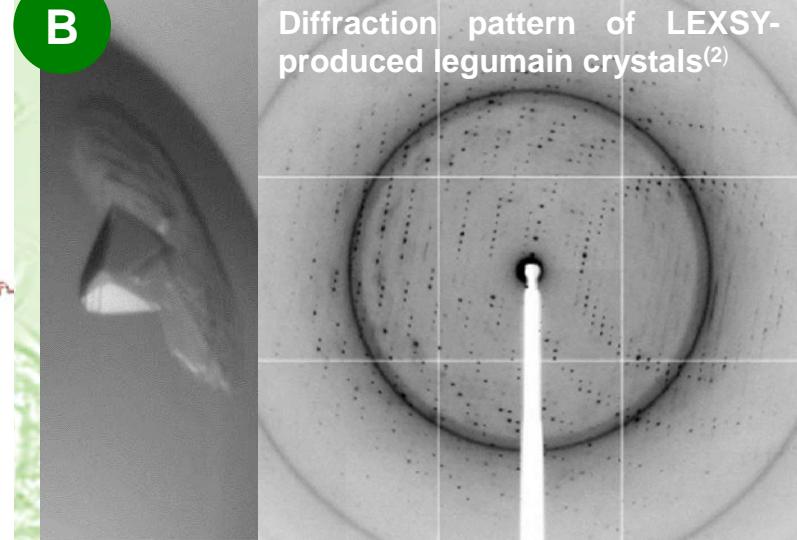
- Crystal structure of LEXSY-made antioxidant human Cu/Zn superoxide dismutase SOD1⁽¹⁾ (A)
- LEXSY-made legumain (cyst. protease) crystals diffract to 2.5 Å⁽²⁾ (B)
- ¹⁵N-HSQC NMR yields assignment of 18x ¹⁵N-Val in a 28 kDa protein⁽³⁾

A



Structure determination of the new P2₁2₁2₁ crystal form of LEXSY-produced human SOD1⁽¹⁾

B



(1) Gazdag et al. (2010) *Acta Crystallographica F*66: 871

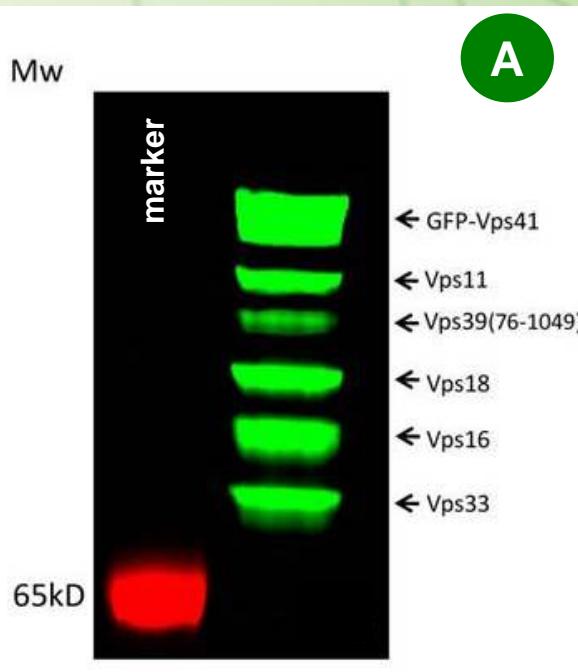
(2) Dall & Brandstetter (2012) *Acta Crystallographica F*68: 24

(3) Niculae et al. (2006) *Protein Expression and Purification* 48: 167

In vitro LEXSY (cell-free expression based on LEXSY cell extracts) is an alternative to traditional wheat germ or RRL systems

In vitro
LEXSY

- LEXSY cell-free production of the to date largest multisubunit membrane protein complex⁽¹⁾ (**A**)
- Fluorescence-based protein interaction studies without purification^(2,3,4)
- Fast high throughput screening^(5,6,7)



Active HOPS membrane tethering complex (**600 kDa**) reconstituted *in vitro* from all six subunits co-expressed by *In vitro* LEXSY⁽¹⁾.

- (1) Guo et al. (2013) PLoS ONE 8: e8153
- (2) Brooks et al. (2014) Science 344: 1249783
- (3) Gambin et al. (2014) eLife 3: e01434
- (4) Han et al. (2014) Journal of Biological Chemistry 289: 7764
- (5) Mureev et al. (2009) Nature Biotechnology 27: 747
- (6) Kovtun et al. (2010) PLOS one 5: e14388
- (7) Kovtun et al. (2011) Methods 55: 58

LEXSY-expressed pharmaceutically relevant enzymes and diagnostics tools

Pathogen
diag-
nostics

- LEXSY-made surface antigens SAG1 and SAG2 of *Toxoplasma gondii* for ELISA kits⁽¹⁾
- LEXSY-made PTX toxin for diagnostics of fungal-caused wheat disease⁽²⁾

Antibodies

- Expression and purification of IgG based scFc fusions⁽³⁾ from LEXSY
- LEXSY-expression and characterization of scFv collection⁽⁴⁾

Pharma
research

LEXSY-made

- Proprotein convertase PC4 (subtilisin kexin) for inhibitor design⁽⁵⁾
- Human liver serine protease Coagulation Factor VII⁽⁶⁾
- Modified human Tissue-Plasminogen Activator (t-PA) that shows >100x higher activity than *E. coli* t-PA⁽⁷⁾
- Glycosylated Amyloid Precursor Protein sAPP α involved in Alzheimer disease⁽⁸⁾

(1) Ebert et al., not published

(2) JBS, not published

(3) Jørgensen et al. (2014) *Microbial Cell Factories* 13: 9

(4) Klatt et al. (2012) *Microbial Cell Factories* 11: 97

(5) Basak et al. (2008) *Protein Expression and Purification* 60: 117

(6) Mirzaahmadi et al. (2011) *Journal of Biomedicine Biotechnology* 2011: 873874

(7) Nazari et al. (2011) *Biotechnology Letters* 33: 503

(8) Klatt et al. (2013) *Journal of Proteome Research* 12: 396

Our LEXSY group...



**Dr. Reinhard
Breitling**
„The LEXSYest Man
Alive“



**Dr. Andreas
Licht**



**Dr. Larissa
Consani
Textor**

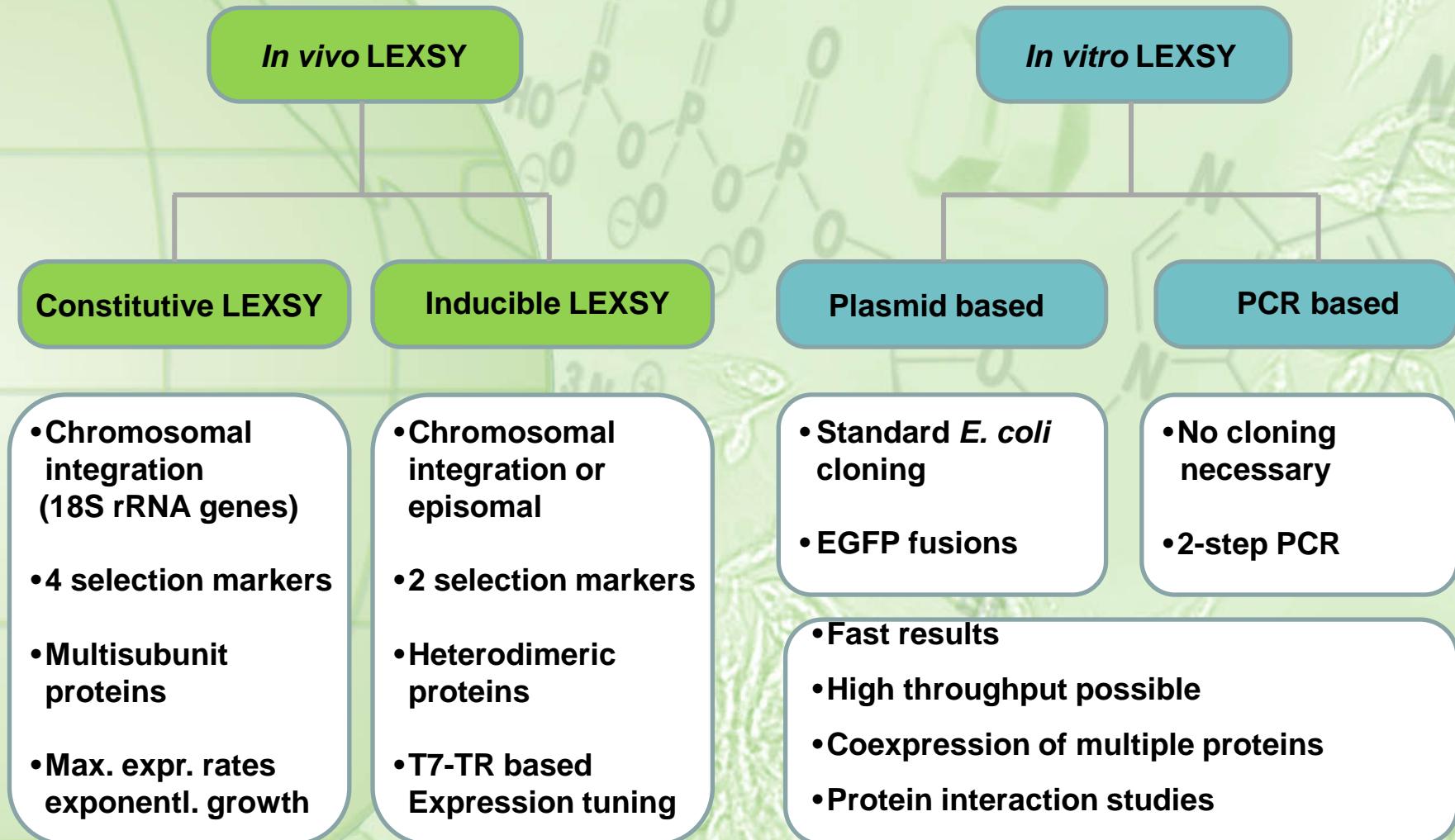


**Stefan
Heiderich**

...answers any question when contacted at expression@jenabioscience.com

backups

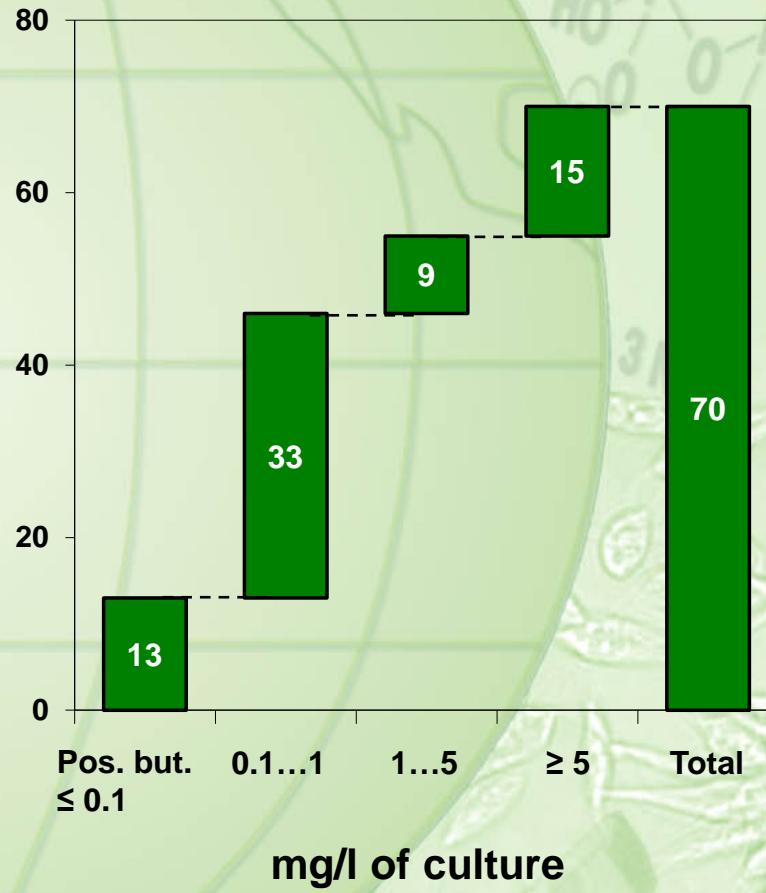
Flexibility of LEXSY expression configurations provides solutions for your needs



80% successful (> 0.1 mg/l) expression projects

Random selection of 70 targets that gave unsatisfactory results in other expression systems

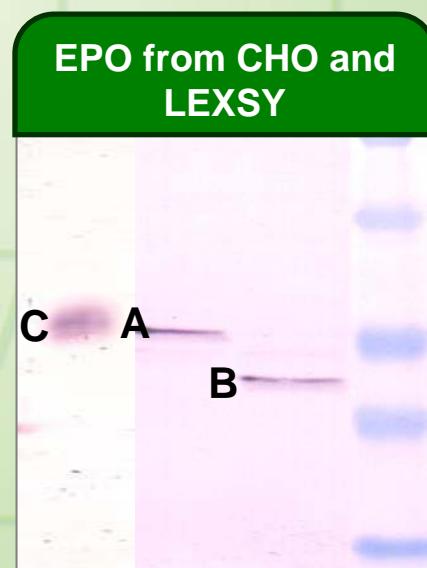
of expression projects



Target protein	Size kDa	Yield mg/L
Cytoplasmic 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

Homogeneously glycosylated human Erythropoietin from LEXSY

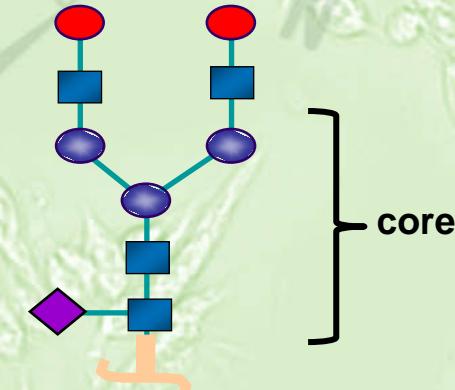
- Completely secreted to the culture medium
- Natively processed at the N-terminus
- Biologically fully active
- Exceptionally homogeneously & mammalian-type N-glycosylated
(biantennary fully galactosylated $\text{Man}_3\text{GlcNAc}_2\text{core}-\alpha-1,6\text{-fucosylated structure}$)



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

Full repertoire of eukaryotic PTMs in LEXSY

PTM	Homol. Targets <i>Leishmania</i>	Heterol. Targets LEXSY	Selected references
Disulfide bond formation (folding)	PMGP46	t-PA	Kahl <i>et al.</i> 1987, Hong <i>et al.</i> 2008 , Soleimani <i>et al.</i> 2007,
Multimerization	RNAP II, RNAP III	Laminin-322 SOD1, TETR	Phan <i>et al.</i> 2009, Gazdag <i>et al.</i> 2010, Kushnir <i>et al.</i> 2005, Martínez-Calvillo <i>et al.</i> 2007
Proteolytic processing -Signal peptide cleavage - Pro-region processing - Furin like processing	Gp63, 3'NT/NU CPB2.8	EPO, IFN γ , SAG1 rPC4 sortilin	Breitling <i>et al.</i> 2002 and unpublished, Brooks <i>et al.</i> 2000, Basak <i>et al.</i> 2008 P. Madsen, not published
Glycosylation - N-Glycosylation - O-Glycosylation - Proteoglycan formation	gp63, sAP, 3'NT/NU sAP, KMP11 sAP	EPO, IFN γ , SAG1	Breitling <i>et al.</i> 2002 and unpublished, Debrabant <i>et al.</i> 2000, Ellis <i>et al.</i> 2002, Ilg <i>et al.</i> 1994, Lippert <i>et al.</i> 1999, McGwire <i>et al.</i> 1996, Olafson <i>et al.</i> 1990, Stierhoff <i>et al.</i> 1998 Weise <i>et al.</i> 2000
Glypiation (GPI anchoring)	Gp63 Proteophosphoglycans MBAP ^{GPI}	rPC4 SAG1	McGwire <i>et al.</i> 1996, Ellis <i>et al.</i> 2002, Foth <i>et al.</i> 2002, Weise <i>et al.</i> (2000) , Basak <i>et al.</i> 2008, Breitling <i>et al.</i> unpublished,
Prenylation	LmLRAB, RAS-CVIM 7 non-assigned 14-140 kDa	Rab7	Chenik <i>et al.</i> 2006, Gillespie <i>et al.</i> 2007, Hasne <i>et al.</i> 1999, Alexandrov <i>et al.</i> not publ.

ARL-1 small G protein ADP-Ribosylation Factor-Like protein 1 of *L. donovani*, **CPB2.8** Cysteine proteinase of *L. mexicana*, **EPO** hu Erythropoietin, **Gp63** most abundant surface glycoprotein, GPI anchored and N-glycosylated HExxH Zn metalloproteinase (leishmanolysin), **H2A** Histone 2A of *L. donovani*, **HSP83-3** Heat shock protein of *L. donovani*, **HSP60** Heat shock protein of *L. donovani*, **IF3** Translation initiation factor 3 subunit of *L. donovani*, **IFN γ** hu intereron gamma, **KMP11** Kinetoplastid membrane protein of *L. donovani*, **LIP2** 60S acidic ribosomal protein P2 of *L. donovani*, **LmLRAB** RAB GTPase of *L. major*, **Lt1200** 1200 kDa Cytoskeletal Giant Protein of *L. tarentolae*, **MBAP^{GPI}** artificially GPI anchored acid phosphatase of *L. mexicana*, **3'NT/NU** Surface Membrane 3'-Nucleotidase/ Nuclease of *L. donovani*, **OADC** bacterial oxaloacetate decarboxylase in *L. mexicana amazonensis*, **PMGP46** promastigote membrane glycoprotein of *L. mexicana amazonensis*, **Rab7** geranylgeranyl transferase component A of *Rattus norvegicus*, **Rbp16** RNA-binding protein of *L. donovani*, **RNAP II** RNA polymerase II of *L.major*, **RNAP III** RNA polymerase III of *L.major*, **rPC4** rat proprotein convertase 4, **S10 & S18** 40S ribosomal proteins of *L. donovani*, **SAG1** surface antigen 1 of *Toxoplasma gondii*, **sAP** secreted acid phosphatase (N- and O-glycosylated and phosphoglycosylation) of *L. mexicana* and *L. donovani*, **SOD1** hu superoxide dismutase, **TETR** tetracycline repressor, **t-PA** hu tissue plasminogen activator, **VG7A5** amastigote spec. Protein of *L. mexicana mexicana*

Full repertoire of eukaryotic PTMs in LEXSY

PTM	Homol. Targets <i>Leishmania</i>	Heterol. Targets LEXSY	Selected references
Methylation	α -Tubulin (D329) Carboxypeptidase (E53) Rbp16 (R104), H2A (S97) ...		Rosenzweig <i>et al.</i> 2008, Jardim <i>et al.</i> 1995
Phosphorylation	Lt1200 (S,T), HSP83-3 (T) RNA helicase II (S) ...		Baqi <i>et al.</i> 2000, Rosenzweig <i>et al.</i> 2008
Acetylation - N-terminal - internal	LIP2 (M1), KMP11 (A2), S10 (S2), IF3 (T2) ... S18 (S15) , HSP60 (E293) ...	SOD1 (A2)	Rosenzweig <i>et al.</i> 2008, Gazdag <i>et al.</i> 2010, Thomas <i>et al.</i> 2009
Myristoylation	ARL-1		Sahin <i>et al.</i> 2008
Amidation	VG7A5		Liu <i>et al.</i> 2000
Glutathionylation	Tb mono-Cys-glutaredoxin 1		Melchers <i>et al.</i> 2007
ADP-ribosylation	TGN-Lysosome trafficking		Sturm <i>et al.</i> 1998 , Sahin <i>et al.</i> 2008
Biotinylation (rec. strain)	-	OADC AVI-EGFP	Detke <i>et al.</i> 2007 Konthur <i>et al.</i> 2009

ARL-1 small G protein ADP-Ribosylation Factor-Like protein 1 of *L. donovani* , **CPB2.8** Cysteine proteinase of *L. mexicana*, **EPO** hu Erythropoietin, **Gp63** most abundant surface glycoprotein, GPI anchored and N-glycosylated HExxH Zn metalloproteinase (leishmanolysin), **H2A** Histone 2A of *L. donovani*, **HSP83-3** Heat shock protein of *L. donovani*, **HSP60** Heat shock protein of *L. donovani*, **IF3** Translation initiation factor 3 subunit of *L. donovani*, **IFN γ** hu intereron gamma, **KMP11** Kinetoplastid membrane protein of *L. donovani*, **LIP2** 60S acidic ribosomal protein P2 of *L. donovani*, **LmLRAB** RAB GTPase of *L. major*, **Lt1200** 1200 kDa Cytoskeletal Giant Protein of *L. tarentolae*, **MBAP^{GPI}** artificially GPI anchored acid phosphatase of *L. mexicana*, **3'NT/NU** Surface Membrane 3'-Nucleotidase/Nuclease of *L. donovani*, **OADC** bacterial oxaloacetate decarboxylase in *L. mexicana amazonensis*, **PMGP46** promastigote membrane glycoprotein of *L. mexicana amazonensis*, **Rab7** geranylgeranyl transferase component A of *Rattus norvegicus*, **Rbp16** RNA-binding protein of *L. donovani*, **RNAP II** RNA polymerase II of *L. major*, **RNAP III** RNA polymerase III of *L. major*, **rPC4** rat proprotein convertase 4, **S10 & S18** 40S ribosomal proteins of *L. donovani*, **SAG1** surface antigen 1 of *Toxoplasma gondii*, **sAP** secreted acid phosphatase (N- and O-glycosylated and phosphoglycosylation) of *L. mexicana* and *L. donovani*, **SOD1** hu superoxide dismutase, **TETR** tetracycline repressor, **t-PA** hu tissue plasminogen activator, **VG7A5** amastigote spec. Protein of *L. mexicana mexicana*



unused

References

- Baqui *et al.* (2000) In Vivo and In Vitro Phosphorylation and Subcellular Localization of Trypanosomatid Cytoskeletal Giant Proteins. *Cell Motility and the Cytoskeleton* **47:** 25
- Basak *et al.* (2008) Recombinant proprotein convertase 4 (PC4) from *Leishmania tarentolae* expression system: Purification, biochemical study and inhibitor design. *Protein Expression and Purification* **60:** 117
- Breitling *et al.* (2002) Non-pathogenic trypanosomatid protozoa as a platform for protein research and production. *Protein Expression and Purification* **25:** 209
- Brooks *et al.* (2000) Processing and trafficking of cysteine proteases in *Leishmania mexicana*. *Journal of Cell Science* **113:** 4035
- Chenik *et al.* (2006) Identification of a new developmentally regulated Leishmania major large RAB GTPase. *Biochem Biophys Res Commun.* **341:** 541
- Debrabant *et al.* (2000) Dissection of the Functional Domains of the *Leishmania* Surface Membrane 3'-Nucleotidase/Nuclease, a Unique Member of the Class I Nuclease Family. *Journal of Biological Chemistry* **275:** 16366
- Detke *et al.* (2007) *Leishmania mexicana amazonensis*: development of a peptide tag useful for labeling and purifying biotinylated recombinant proteins. *Exp Parasitol.* **115:** 221
- Ellis *et al.* (2002) Processing and Trafficking of *Leishmania mexicana* GP63. *Journal of Biological Chemistry* **277:** 27968
- Foth *et al.* 2002 *Leishmania* major proteophosphoglycans exist as membrane-bound and soluble forms and localise to the cell membrane, the flagellar pocket and the lysosome. *Int J Parasitol.* **32:**1701
- Gazdag *et al.* (2010) Purification and crystallization of human Cu/Zn superoxide dismutase recombinantly produced in the protozoan *Leishmania tarentolae*. *Acta Crystallographica F* **66:** 871
- Gillespie *et al.* (2007) C-terminal proteolysis of prenylated proteins in trypanosomatids and RNA interference of enzymes required for the post-translational processing pathway of farnesylated proteins. *Mol Biochem Parasitol.* **2007 153:** 115
- Hasne *et al.* (1999) Characterization of prenylated protein methyltransferase in *Leishmania*. *Biochem. J.* **342:** 513
- Hong *et al.* (2008) Identification and enzymatic activities of four protein disulfide isomerase (PDI) isoforms of *Leishmania amazonensis*. *Parasitol. Res.* **102:** 437

References

- Ilg *et al.* (1994) O- and N-Glycosylation of the *Leishmania mexicana*-secreted Acid Phosphatase. *Journal of Biological Chemistry* **269:** 24073
- Jardim *et al.* (1995) Isolation and structural characterization of the *Leishmania donovani* kinetoplastid membrane protein-11, a major immunoreactive membrane Glycoprotein. *Biochem. J.* **305:** 307
- Kahl *et al.* (1987) Structural and antigenic characterization of a species- and promastigote-specific *Leishmania mexicana amazonensis* membrane protein. *J. Immunol.* **138:**1587
- Konthur *et al.* 2009 *In Vivo* Biotinylation of Recombinant Proteins in the Protozoan Host *Leishmania tarentolae*. PEPTalk 2009 January 11 - 16, Coronado, CA
- Kushnir *et al.* (2005) Development of an inducible protein expression system based on the protozoan host *Leishmania tarentolae*. *Protein Expression and Purification* **42:** 37
- Lippert *et al.*(1999) Phosphoglycosylation of a secreted acid phosphatase from *Leishmania donovani*. *Glycobiology*. **9:** 627
- Liu *et al.* (2000) Isolation and analysis of a new developmentally regulated gene from amastigotes of *Leishmania mexicana mexicana*. *Parasitol Res.* **86:** 140
- Martínez-Calvillo *et al.* (2007) Characterization of the RNA polymerase II and III complexes in *Leishmania major*. *Int J Parasitol.* **37:** 491
- McGwire *et al.* (1996) Posttranslational Regulation of a *Leishmania* HEXXH Metalloprotease (gp63). *Journal of Biological Chemistry* **265:** 12240
- Melchers *et al.* 2007 Glutathionylation of trypanosomal thiol redox proteins. *J Biol Chem.* **282:** 8678
- Olafson *et al.* (1990) Structures of the N-linked oligosaccharides of Gp63, the major surface glycoprotein, from *Leishmania mexicana amazonensis*. *Journal of Biological Chemistry* **269:** 24073
- Phan *et al.* (2009) The production of recombinant human laminin-332 in a *Leishmania tarentolae* expression system. *Protein Expression and Purification* **68:** 79
- Rosenzweig *et al.* (2008) Post-translational modification of cellular proteins during *Leishmania donovani* differentiation. *Proteomics*. **8:** 1843
- Sahin *et al.* (2008)The *Leishmania* ARL-1 and Golgi Traffic. *PLoS One.* **3:** e1620

References

- Soleimani *et al.* (2007) Expression of human tissue-type plasminogen activator (t-PA) in *Leishmania tarentolae*. *Biotechnology & Applied Biochemistry* **48**: 55
- Sturm *et al.* (1998) Characterization of a GTP-binding protein in the ADP-ribosylation factor subfamily from *Leishmania tarentolae*. *Biochim Biophys Acta* **1442**: 347
- Stierhoff *et al.* (1998) Structure of a filamentous phosphoglycoprotein polymer: the secreted acid phosphatase of *Leishmania mexicana*. *J. Mol. Biol.* **282**: 137
- Thomas *et al.* (2009) Histone acetylations mark origins of polycistronic transcription in *Leishmania major*. *BMC Genomics* **10**: 152
- Weise *et al.* (2000) Distribution of GPI-anchored proteins in the protozoan parasite *Leishmania*, based on an improved ultrastructural description using high-pressure frozen cells. *Journal of Cell Science* **113**: 4587