

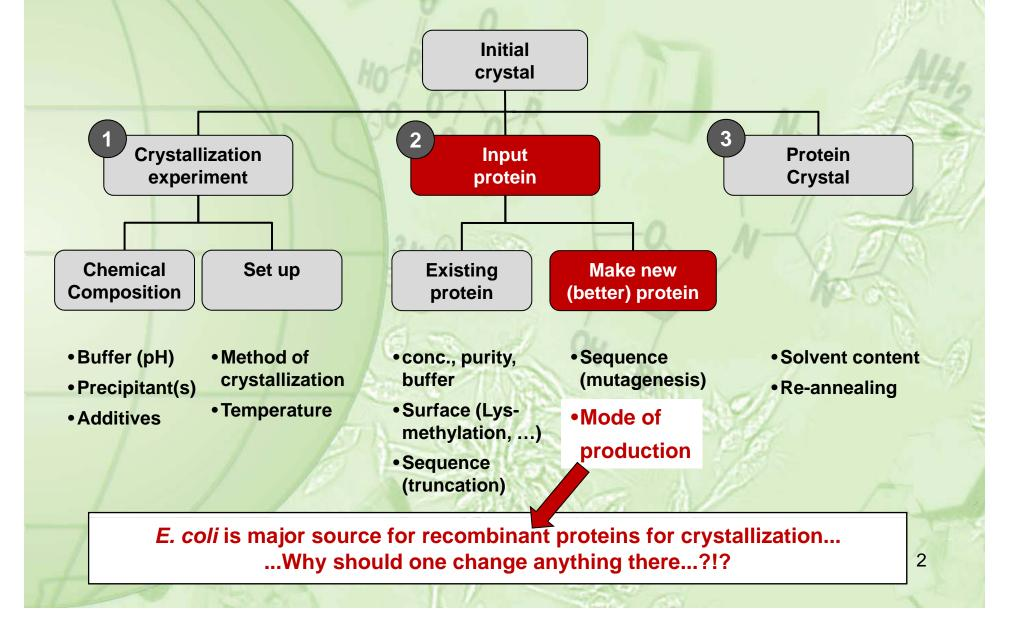


# From gene to crystallization within two days

**LEXSY** cell-free protein production

Jena Bioscience GmbH Löbstedter Str. 80 07749 Jena, Germany Tel.: +49-3641-628-5000 Fax: +49-3641-628-5100 e-Mail: info@jenabioscience.com

## There are more options for optimising hits than adjusting pH Protein sample equally important: No good protein → no good crystal



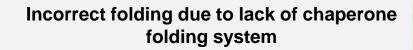
## Two major reasons why *E. coli* proteins often do not crystallize *E. coli* in general not a good host for expression of "interesting" proteins

Typical "symptoms" with *E. coli* derived

## proteins

### Underlying shortcoming of *E. coli*

- Insoluble / Inclusion bodies
- Aggregate-forming → precipitates over time and compromised biological activity
  - Low solubility, e.g. cannot be concentrated (< 10 mg/ml)</li>
    - Unstable/fast degradation



→ Applies to almost all proteins<sup>(1)</sup> of higher organisms expressed in *E. coli* 

- Low expression yields (< 1 mg/L)
  - Unstable/fast degradation
- 2
- Compromised biological activity and/or ligand binding
  - Low solubility

Incorrect processing due to lack of posttranslational modification<sup>(2)</sup> machinery

→ Applies to almost all proteins<sup>(1)</sup> of higher organisms expressed in *E. coli* 

3

What should one do when low quality protein is obtained from E. coli...?

(1) >50% of human proteins are glycosylated, 10<sup>6</sup> protein variants present in humans (Walsh & Jefferis, 2006, Nat. Biotech. 24: 1241
 (2) PTMs: glycosylation<sup>(1)</sup>, disulfide bond formation, phosphorylation, acetylation, sulfation, amidation, prenylation, ...)

## Changing mode of protein production – a difficult trade-off But often the better choice than "riding a dead horse<sup>(1)</sup>"

Being faithful to *E. coli* 

Changing mode of production

Plenty of opportunities to optimize crystallization experiment, i.e. riding the dead horse<sup>(1)</sup> Starting again from scratch

- (Sub)cloning gene of interest
- Creation of expression strain
- Preparative cultivation/expression
- Protein harvest & purification

... or something completely different...

Changing mode of production may sometimes be the wiser strategy...

...and not as cumbersome as thought

(1) Dakota tribal wisdom, see http://www.tysknews.com/LiteStuff/riding a dead horse.htm

## May I introduce LEXSY: A Combination of robustness and eukaryotic features





## Easy construction and rapid growth of LEXSY expression strains

- Biosafety level 1 (as E. coli)
- Plasmid generation in *E. coli* shuttle vectors
- Cultivation in inexpensive media at 26°C
- Expression yields of up to 300 mg/l of culture

## Eukaryotic protein folding and modification machinery

- Chaperones for folding (no inclusion bodies)
- Full range of PTMs (including exceptionally homogenous mammaliantype N-glycosylation)

### Constitutive or inducible, intracellular or secretory versions

Coexpression of up to 4 target genes

Does it work for X-ray crystallography...?

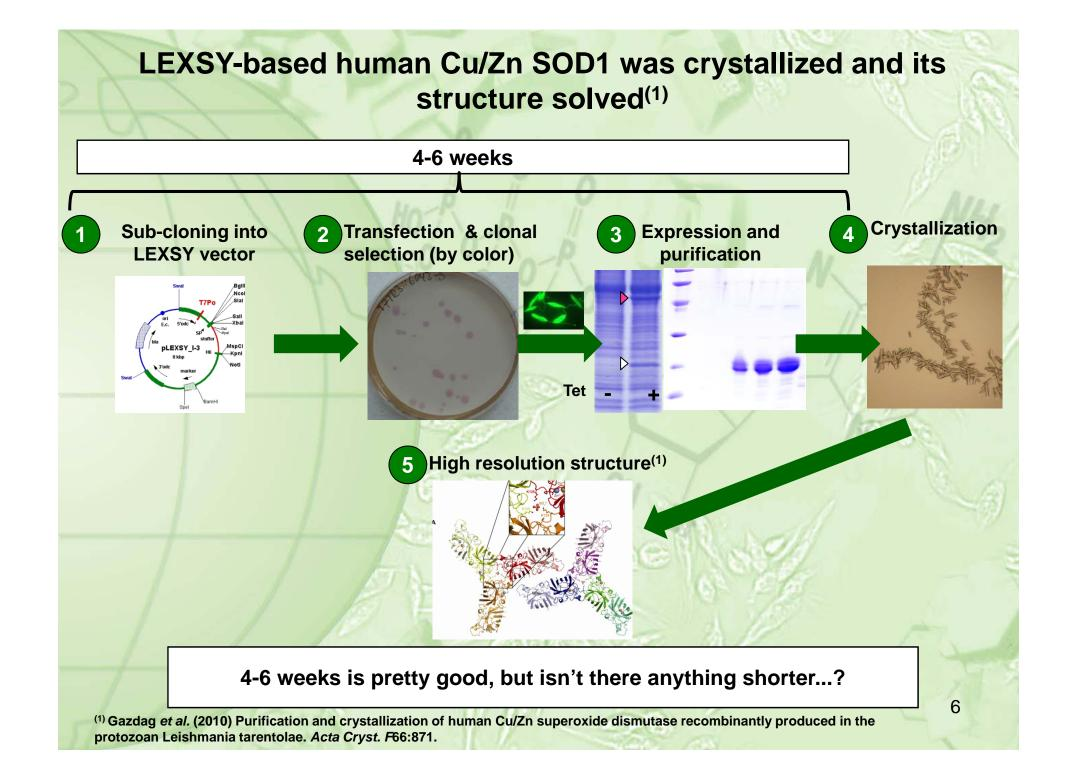
#### EPO from:

CHO LEXSY (smear) (single species)

(1)

(1) Deglycosylated

5

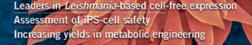


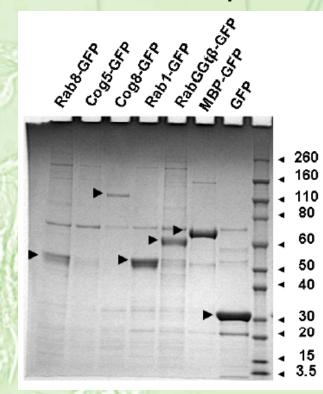
## Yes there is: LEXSY in vitro expression kit now available



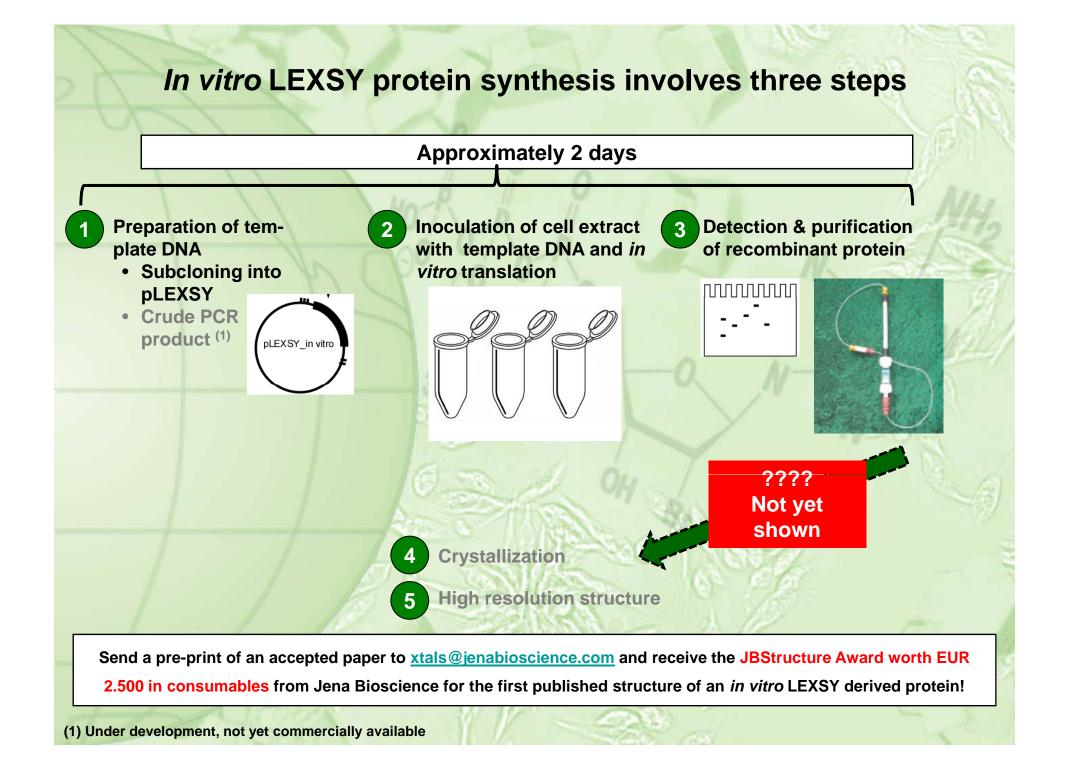
- Based on LEXSY transcription-translationcompetent cell extracts
- No cell culturing necessary

Yields of 0.3 mg/ml of extract within 2 hours after initiation of *in vitro* expression





7



## Summary: Two straightforward choices for skipping E. coli

#### In vivo expression



Programmable LEXSY (S1) cells (here: eGFP)

#### Yields up to 300 mg/L of culture

Proven for structure determination (SOD1, eGFP)

#### Advantages:

- Eliminates all shortcomings of E. coli
- Easy to handle
- No cell culture equipment necessary

#### **Disadvantage:**

 Requires going back to wet lab (cloning → transfection & cultivation → protein purification, 4-6 weeks total)

#### In vitro expression (cell-free)



Programmable cell-extracts (*in vitro*)

Yields up to 0.3 mg/ml of cell extract

Not yet proven for structure determination

#### Advantages:

- Eliminates all shortcomings of E. coli
- Easy to handle
- Very fast results

#### **Disadvantage:**

- Lower yields (0.3 mg/ml of cell extract ) and more expensive consumables
- → But 0.1 mg of protein allows crystal set ups in roughly 60 wells<sup>(1)</sup>

(1) Assuming a protein solution of 10 mg/ml and 150 nl per well

## Acknowledgements



Kirill Alexandrov & group

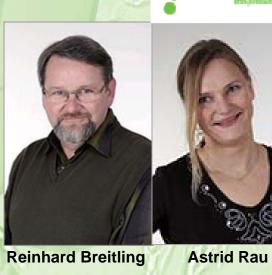


Wulf Blankenfeld & group

Jana Blosciance



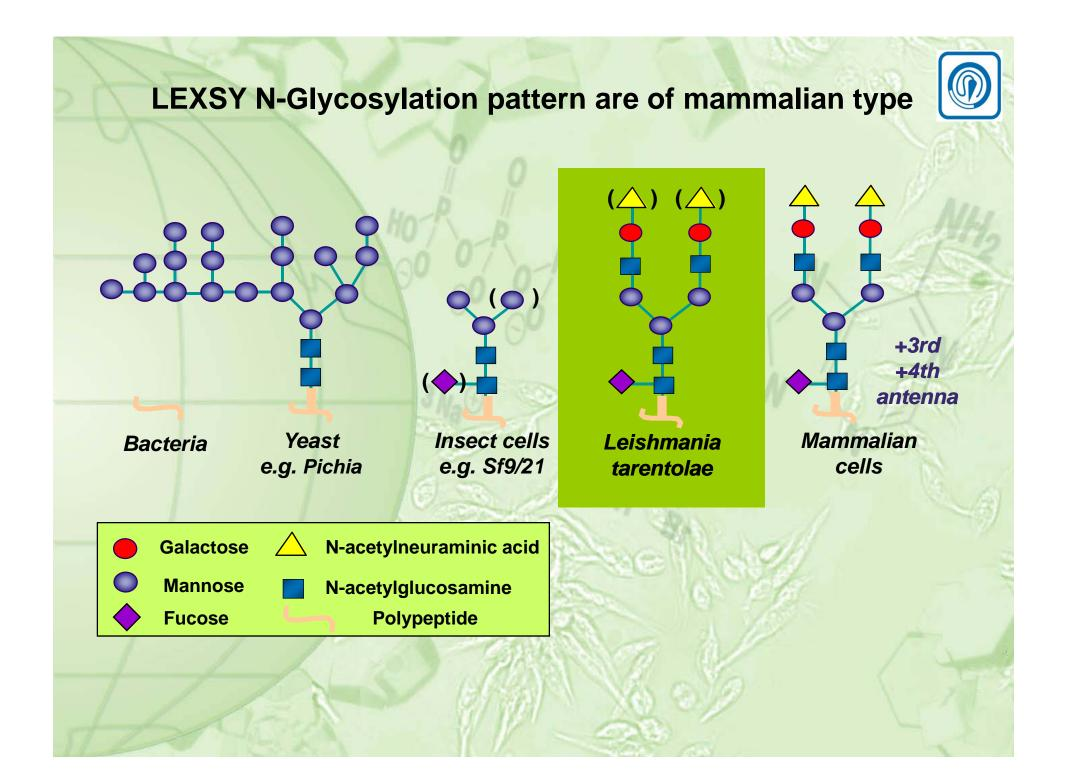
Peter Bayer



**Thomas Billert** 

## unused

11



## "JBStructure Award": EUR 2.500 in consumables for first published structure of *in vitro* LEXSY protein

#### In vitro expression (cell-free)



Programmable cell-extracts (*in vitro*)

Yields up to 0.3 mg/ml of cell extract

Not yet proven for structure determination

Send a pre-print of an accepted paper to <u>xtals@jenabioscience.com</u> and receive the JBStructure Award worth EUR 2.500 in consumables from Jena Bioscience for the first published structure of an *in vitro* LEXSY derived protein!

## Again – commercially available cell-free systems have shortcomings



- Lack of chaperone based mechanisms
- Absence of posttranslational modifications





Rabbit

reticulocyte

eato

14

Difficult scale up

- Laborious and expensive preparation
  - Difficult to manipulate genetically

## Shortcomings of conventional expression systems require an alternative expression system



- Insufficient folding of complex proteins of higher organisms
- Lack of posttranslational modifications
- Endotoxins



- 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



- Long
  developmental
  cycles
- Complex
  downstream
  processing
- Contamination
  problems

Current expression systems are not ready for proteomics era Need eukaryotic machinery but prokaryotic robustness