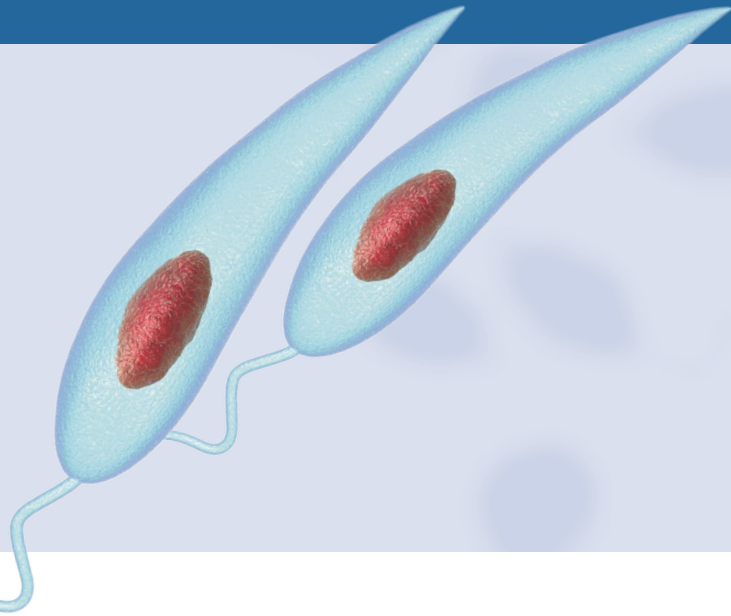


LEXSY

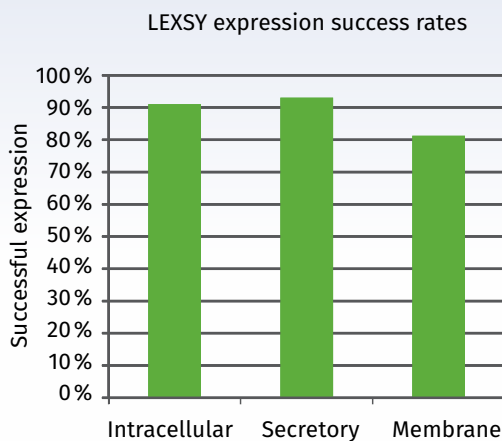
Eukaryotic Protein Expression



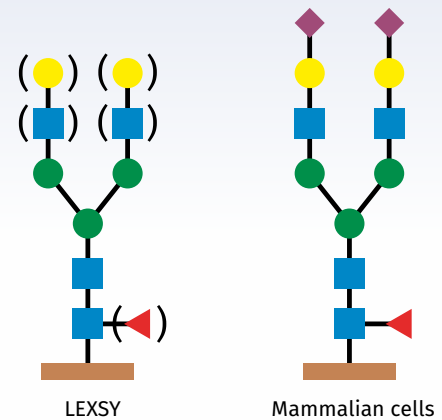
LEXSY is a eukaryotic protein expression system based on the S1-classified unicellular organism *Leishmania tarentolae*. It features...

- › Simple Handling
- › Eukaryotic Posttranslational modifications
- › High success rates for difficult proteins
- › An S1-classified expression host

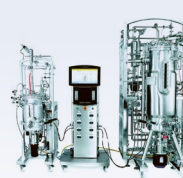
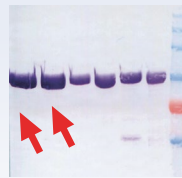
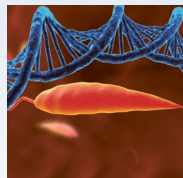
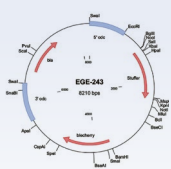
LEXSY achieves high success rates, even with difficult-to-express proteins



LEXSY performs mammalian-type glycosylation



LEXSY is fast: From initial cloning to purified protein in 6 weeks



Expression plasmid construction

1 week

Transfection & selection

2 weeks

Expression evaluation

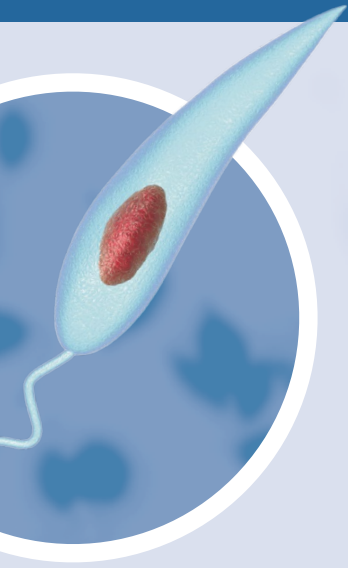
1 week

Scale-up

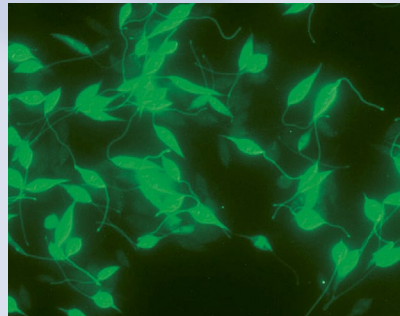
1 week

Protein purification

1 week



LEXSY performs a wide range of eukaryotic postranslational modifications, including:



Leishmania cells expressing enhanced GFP

- › N- and O-glycosylation
- › Disulfide bond formation
- › Localization / Secretion signal processing
- › Proteolytic processing
- › Phosphorylation
- › Lipidation
- › Acetylation

LEXSY starter kits supply you with everything you need for your first expression run

Expression vector

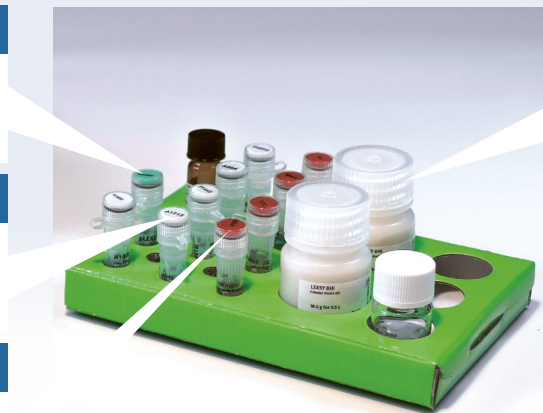
For intracellular or secretory expression, four markers available

LEXSY primers

Sequencing and diagnostic primer pairs

LEXSY antibiotic

For selection of recombinant strains



LEXSY medium

All components for 1 litre of growth medium

LEXSY host

Glycerol stocks



Custom expression service:

Jena Bioscience offers expression of your protein using our proprietary expression system LEXSY.

We offer:

- › High success rates
- › Speed and flexibility
- › Modular work plans - tailored to your needs
- › Attractive cost schemes

