

LEXSInduce3 Expression Kit (contains the integrative vector

pLEXSY_I-neo3)

The LEXSInduce3 kit contains one of three expression vectors of choice which differ in the antibiotic resistance marker and the option to monitor expression online by a fluorescence reporter. The vector pLEXSY_I-neo3 (encoding aminoglycoside phosphotransferase) allows selection with the antibiotic LEXSY Neo, whereas pLEXSY_I-ble3 (encoding bleomycin resistance gene) permits selection with the antibiotic LEXSY Bleo. The vector pLEXSY_I-blecherry3 facilitates selection with the antibiotic LEXSY Bleo and, in addition, offers the possibility to monitor induction during cultivation by coexpressed fluorescence. The control vector provided in the kit contains the EGFP gene linked to the respective selection marker.

The pLEXSY_I-3 vectors enable inducible expression of target proteins either with or without secretory signal peptide provided on the vectors. For customer convenience the same vector can be used for cloning of ORFs either for cytosolic or for secretory expression. The LmSAP secretory signal peptide encoded on these vectors was derived from the gene for secreted acid phosphatase (Lmsap1) of *Leishmania mexicana*. In-frame fusion of the ORF of a target protein to this signal peptide allows secretory expression in LEXSY hosts, whereas cloning into any of the restriction sites at the 5' end of the signal peptide-encoding sequence will result in cytosolic expression. An illustration of the inducible LEXSY can be downloaded from the Jena Bioscience website at

http://www.jenabioscience.com/images/b3e879b381/Illustration_inducible_LEXSY.pdf

One main advantage of LEXSY is the mammalian-type posttranslational modification of target proteins, such as glycosylation, phosphorylation or prenylation. Recombinant human erythropoietin (EPO) purified at Jena Bioscience from LEXSY was biologically active, natively processed at the N-terminus, and N-glycosylated. The N-glycosylation profile was exceptionally homogeneous, with a biantennary oligosaccharide and the Man3GlcNAc2 core structure accounting for >90% of the glycans present. *L. tarentolae* is thus the first described biotechnologically useful unicellular eukaryotic host producing biantennary, fully galactosylated, core- α -1,6-fucosylated N-glycans. This N-glycosylation profile was coincident with the profile of recombinant human Interferon- γ expressed in LEXSY and of LEXSY host Gp63 glycoprotein. An illustration of LEXSY glycosylation can be downloaded from the Jena Bioscience website at

http://www.jenabioscience.com/images/b3e879b381/Illustration_LEXSY_Glycosylation.pdf

Ordering Information

[Order Online »](#)

JBS-EGE-1410neo	1Kit
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Contents

- LEXSY host T7-TR: kit contains three vials with 1.6 ml each of frozen glycerol stocks of LEXSY host T7-TR.
- pLEXSY_I-3 expression vector pLEXSY_I-neo3 (Cat. No. JBS-EGE-245), 5 µg in 10 mM Tris HCl pH 8.0
- pLEXSY_I-3 control plasmid with EGFP gene pLEXSY_I-egfp-neo3 (Cat. No. JBS-EGE-248), 5 µg in 10 mM TrisHCl pH 8.0
- Primers for diagnostic PCR and sequencing:
neo forward primer A1432, for pLEXSY_I-neo3
Insert sequencing reverse primer A264
5`utr (aprt) reverse primer A1715
odc forward primer A1304
odc reverse primer P1510
- LEXSY BHI, Powder for preparation of liquid cultivation medium
- LEXSY Neo* for pLEXSY_I-neo3, 1 ml ready-to-use 1000x stock solution, 50 mg/ml, filter-sterilized
- inducer of the T7-TR system: LEXSY Tet (Tetracycline), 1 ml ready-to-use 1000x stock solution, 10 mg/ml, filter sterilized
- Hemin, 2 ml ready-to-use 500x stock solution, steril-filtered
- Pen-Strep, 5 ml ready-to-use 200x stock solution, steril-filtered

Technical Info / Product Notes

For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com