

## LEXSY Cultivation Kit T7-TR

*contains LEXSY host T7-TR*  
for inducible LEXSY expression vectors

Cat.-No.	Amount
LT-111	1 Kit

### Description:

The **LEXSY Cultivation Kit T7-TR** was developed for initial cultivation and establishment of *Leishmania tarentolae* strain T7-TR in suspension culture. The kit is shipped on dry ice.

### The kit contains:

3 vials with 1.6 ml each of frozen glycerol stocks of LEXSY host T7-TR (*Leishmania tarentolae* strain T7-TR expressing bacteriophage T7 RNA polymerase and TET repressor)

50 ml LEXSY BHI liquid medium complete

For initial inoculation; contains Hemin, LEXSY NTC, LEXSY Hygro and PenStrep

37 g LEXSY BHI powder

2x 18.5 g for preparation of 2x 500 ml LEXSY BHI growth medium

2 ml Hemin 500x stock solution

0,25% solution of porcine Hemin in 50% Triethanolamine for 1 L LEXSY BHI medium

5 ml Pen-Strep stock solution, 200x

10.000 units of penicillin (base) and 10.000 µg of streptomycin (base)/ml in 0.85% saline; for 1 L LEXSY BHI medium

1 ml LEXSY NTC and 1 ml LEXSY Hygro

for maintenance of T7 polymerase and TET repressor genes.

4 cryo vials with each 0.4 ml 80% glycerol, sterile sterile, for preparation of glycerol stocks

3 culture flasks for 10 ml medium

For initial inoculation.

### Storage conditions:

*Glycerol stocks of LEXSY host*

**Upon arrival the glycerol stocks must be stored at -80°C or inoculated into LEXSY BHI provided! Do not freeze-thaw the stocks!**

*LEXSY BHI liquid medium*

Store at 4°, stable for 2 weeks

*LEXSY BHI powder*

Store at ambient temperature, stable for 3 years

*Hemin stock solution*

Store at 4°C in the dark, stable for 6 months

*LEXSY NTC, LEXSY Hygro and Pen-Strep stock sol.*

Store at -20°C, stable for 6 months

*cryo vials with glycerol*

store at ambient temperature, stable for 6 months.

### Reactivation of LEXSY host:

Thaw glycerol stock on ice and inoculate the **entire content** of the vial into 10 ml of LEXSY BHI medium provided. Incubate at 26°C and dilute as required.

### Preparation of LEXSY BHI growth medium:

Dissolve 37 g/l LEXSY BHI powder in deionized water and autoclave 15 min at 121°C. Add Hemin, LEXSY NTC, LEXSY Hygro and PenStrep. Store at 4°C and use within two weeks.

### Preparation of glycerol stocks:

Add 1.2 ml of growing culture (ca.  $6 \times 10^7$  cells/ml) to one vial with 0.4 ml 80% glycerol, mix, incubate 10 min at RT, 1 h on ice and o/n at -20°C. Transfer to -80°C. Strains can be stored this way for several years.

Find more information in the User's Guide for *Leishmania tarentolae* Gene Expression Starter Kits sect. 3

**FOR RESEARCH USE ONLY.**

**NOT INTENDED FOR HUMAN OR ANIMAL  
DIAGNOSTIC OR THERAPEUTIC USES.**