



LEXSY Cultivation Kit P10

Contains LEXSY host P10
For constitutive expression

Cat. No.	Amount
LT-102	1 Kit

For research use only! Not intended for human or animal diagnostic or therapeutic uses.

Shipping: shipped on dry ice/on gel packs

Storage Conditions: Store components as indicated on individual labels

Shelf Life: See individual components

Description:

The LEXSY Cultivation Kit P10 was developed for initial cultivation and establishment of *Leishmania tarentolae* laboratory strain P10 in suspension culture.

Content:

3 vials with 1.6 ml of frozen glycerol stocks of LEXSY host P10 each (*Leishmania tarentolae* laboratory strain P10)

Upon arrival the glycerol stocks must be stored at -80 °C or inoculated into the provided LEXSY BHI liquid medium. Do not thaw and re-freeze the stocks.

50 ml LEXSY BHI liquid medium complete

For initial inoculation, contains Hemin and PenStrep.

Store at 4 °C, stable for 2 weeks

37 g LEXSY BHI powder

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

Store at ambient temperature, stable for 3 years

2 ml Hemin stock solution, 500x

0,25 % solution of porcine Hemin in 30 % Triethanolamine for 1 L LEXSY BHI medium.

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

5 ml Pen-Strep stock solution, 200x

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

Store at -20 °C, stable for 6 months

4 cryo vials with 0.4 ml 80 % glycerol each, sterile

For preparation of glycerol stocks.

Store at ambient temperature, stable for 6 months

3 cell culture flasks for 10 ml medium

For initial inoculation.

Biosafety level:

1, Non-pathogenic

Source:

Tarentola annularis

Reactivation of LEXSY host:

Thaw glycerol stock on ice for ca. 20 minutes and inoculate the entire content of the vial into 10 ml of LEXSY BHI medium in one of the provided cell culture flasks. Incubate at 26 °C in the dark and dilute as required.

Preparation of LEXSY BHI growth medium:

Dissolve 37 g/l LEXSY BHI powder in deionized water and autoclave max. 15 min at 121 °C. Add Hemin and PenStrep to 1x final concentration. Store at 4 °C in the dark and use within two weeks.



LEXSY Cultivation Kit P10

Contains LEXSY host P10
For constitutive expression

Preparation of glycerol stocks:

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