



## LEXSY BHI - Liquid Media Kit

sterile, brain-heart infusion-based medium  
recommended for transfection and strain maintenance

Cat. No.	Amount
ML-411S	1 L
ML-411L	5 x 1 L

**For general laboratory use.** Not intended for human or animal diagnostic or therapeutic uses.

**Shipping:** Shipped at ambient temperature/on gel packs

**Storage Conditions:** Store components as indicated on individual labels

**Shelf Life:** 12 months

### Applications:

For the growth of LEXSY hosts P10 (Cat.-No. LT-101) and T7-TR (Cat.-No. LT-110).

### Description:

Three-component Kit for preparation of complex *Leishmania* cultivation medium.

### Content:

*LEXSY BHI, liquid*

*Hemin stock solution, 500x* (Cat.-No. ML-108):

0,25 % solution of Hemin in 30 % Triethanolamine, sterile

*Pen-Strep stock solution, 200x* (Cat.-No. ML-105):

10.000 U/ml of penicillin G sodium salt and 10.000 µg/ml of streptomycin sulfate in 0.85 % saline, sterile

Cat. No.	Content	Amount
<b>ML-411S</b>	1 l medium 2 ml Hemin 5 ml Pen-Strep	for 1 l medium
<b>ML-411L</b>	5 x 1 l medium 10 ml Hemin 25 ml Pen-Strep	for 5 l medium

### Storage conditions:

*LEXSY BHI, liquid:*

Store at room temperature

Stable for 12 months

*Hemin stock solution, 500x:*

Store at 4 °C in the dark

Stable for 12 month

*Pen-Strep stock solution, 200x:*

Store at -20 °C

Stable for 12 month

### Preparation of complete *Leishmania* cultivation medium:

Add to 1 l LEXSY BHI:

5.0 ml Pen-Strep solution

2.0 ml Hemin stock solution to a final concentration of 5 µg/ml

Final pH = 7.6 ± 0.2

Store in the dark at 4 °C.

Completed medium is useable for up to 2 weeks. If the completed medium is to be used after this period, appropriate amounts of Hemin and Pen-Strep have to be re-added.



## LEXSY BHI - Liquid Media Kit

sterile, brain-heart infusion-based medium  
recommended for transfection and strain maintenance

### Selected References:

[1] Zauner *et al.* (2018) Structural Analyses of *Arabidopsis thaliana* Legumain  $\gamma$  Reveal Differential Recognition and Processing of Proteolysis and Ligation Substrates. *J Biol Chem* **293**: 8934