



LEXSY BHI - Powder Media Kit

brain-heart infusion-based
recommended for transfection and strain maintenance

Cat. No.	Amount
ML-412S	for 1 L
ML-412L	for 5 L
ML-412XL	for 10 L
ML-412XXL	for 50 L

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

Shipping: shipped 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 complete *Leishmania* cultivation medium.

Content:

LEXSY BHI, powder

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
ML-412S	37 g powder 2 ml Hemin 5 ml Pen-Strep	for 1 l medium
ML-412L	185 g powder 10 ml Hemin 25 ml Pen-Strep	for 5 l medium
ML-412XL	370 g powder 20 ml Hemin 50 ml Pen-Strep	for 10 l medium
ML-412XXL	1850 g powder 5x 20 ml Hemin 250 ml Pen-Strep	for 50 l medium

Storage conditions:

LEXSY BHI, powder:

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:

Dissolve 37 g/l LEXSY BHI powder in deionized water and autoclave for 15 min at 121 °C. Control temperature and time with a reference

**LEXSY BHI - Powder Media Kit**

brain-heart infusion-based
recommended for transfection and strain maintenance

with the same volume, since overexposure leads to de-composition of ingredients and a dark brown color of the medium, which adversely affects growth of LEXSY strains. Proper autoclaved medium is of amber color. Store medium at room temperature before addition of Hemin and Pen-Strep (stable for 12 months). As an alternative, the medium can also be sterile filtered.

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.

Selected References:

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