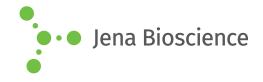
### **DATA SHEET**





#### ■ LEXSY Plating Kit basic

components for solid medium brain-heart infusion-based, without FCS, NC membranes, spatula, dishes & serological pipettes

Cat. No.	Amount
ML-453	for 40 plates

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

Shipping: Shipped at ambient temperature/on gel packs

 $\textbf{Storage Conditions:} \ \ \text{store components as indicated on individual}$ 

labels

Shelf Life: 12 months

#### **Description:**

Core kit for preparation of solidified medium for clonal selection of Leishmania tarentolae. Plasticware and inactivated bovine serum (iFCS) to be provides by customer.

#### Content:

- 2 x 250 ml of Agar base
- 400 ml of buffered liquid medium:
   This component is provided in excess to allow for individual batches.
- 2 ml of Hemin stock solution, 500x Cat.-No. ML-108S:
   0,25% solution of Hemin in 30% Triethanolamine, sterile
- 5 ml of Pen-Strep stock solution, 200x Cat.-No. ML-105S: 10.000 units of penicillin (base) and 10.000 µg of streptomycin (base)/ml as penicillin G sodium salt and streptomycin sulfate in 0.85% saline, sterile

#### All components are provided sterile.

This kit does **not contain** inactivated bovine serum (iFCS), nitrocellulose membranes, spatula, dishes, pipettes which are provided with the **LEXSY Plating Kit comfort** (Cat.-No. ML-451).

#### Storage conditions:

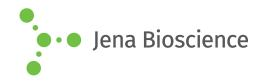
- Agar base:
  - Store at room temperature Stable for 12 months
- Buffered liquid medium:
  - Store at 4 °C
  - Stable for 12 month
- Hemin stock solution, 500x:
  - Store at 4 °C in the dark
  - Stable for 12 month
- Pen-Strep stock solution, 200x: Store at -20 °C
  - Stable for 12 month

## Preparation of 500 ml LEXSY plating medium (20 plates, each 25 ml):

- Microwave-melt 250 ml Agar base in provided glass flask and keep at 55 °C.
- Add pre-warmed (37 °C) buffered liquid medium to ca. 400 ml.
- Add 50 ml iFCS (not provided), 1 ml Hemin stock and 2.5 ml Pen-Strep.
- Fill with buffered liquid medium to 500 ml.
- Add appropriate antibiotics (not provided).
- Pour plates with 25 ml pre-warmed (55 °C) serological pipette, avoid air bubbles.
- After solidification, dry the plates open under the laminar flow



## **DATA SHEET**





# LEXSY Plating Kit basic

components for solid medium brain-heart infusion-based, without FCS, NC membranes, spatula, dishes & serological pipettes

box for exactly 10 min, than store them closed bottom-up.
After plating, seal plates with parafilm and incubate bottom- up
for 5 - 9 days at 26 °C until colonies appear.

Prepare the plates freshly on the day of plating!