

LEXSY Plating Kit comfort

with nitrocellulose membranes, spatula, dishes, pipettes

Cat.-No.	Amount
ML-451	1 liter medium for 40 plates

Description:

All-inclusive kit for preparation of solidified medium for clonal selection of *Leishmania tarentolae*.

The kit contains:

2 x 250 ml Agar base.

400 ml buffered liquid medium.

This component is provided in excess to allow for individual batches.

2 x 50 ml heat-inactivated fetal calf serum (iFCS).

2 ml Hemin stock solution, 500x (ML-108):
0,25% solution of porcine Hemin in 50% Triethanolamine.

5 ml Pen-Strep stock solution, 200x (ML-105):
10.000 units of penicillin (base) and 10.000 µg of streptomycin (base)/ml in 0.85% saline.

40 x Nitrocellulose membranes diameter 82 mm.
For easy plating, fluorescence monitoring and transfer of colonies to alternative conditions (e.g. induction of the LEXSY T7 polymerase-based gene expression).

40 x Petri dishes diameter 90 mm.

40 x Plating spatula in single bags.

Serological pipettes 4 x 25 ml, 2 x 10 ml,
2 x 2 ml.

All components are provided sterile

Storage conditions:

Agar base

Store at room temperature. Stable for 6 months.

Buffered liquid medium

Store at 4°. Stable for 6 months

iFCS

Store < -20°C. Stable for 6 months.

Hemin stock solution

Store at 4°C in the dark. Stable for 6 months.

Pen-Strep stock solution

Store at -20°C. Stable for 6 months.

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

Microwave-melt 250 ml Agar base in provided glass flask keep at 55°C.

Add pre-warmed (37°C) buffered liquid medium to ca. 400 ml.

Add 50 ml iFCS, 1 ml Hemin stock and 5 ml PenStrep.

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 box for exactly 10 min, then store them closed bottom-up.

Prepare the plates freshly on the day of plating!

After plating, seal plates with parafilm and incubate bottom-up for 5-7 days at 26°C until colonies appear.

Find more information in the User's Guide for *Leishmania tarentolae* Gene Expression Starter Kit (download pdf from the LEXSY section of our website).

FOR RESEARCH USE ONLY.

Not intended for human or animal diagnostic or therapeutic uses.