

LEXSY YS - Powder Media Kit

yeast-soybean based, casein-free, for medium free of animal components
 (incl. Glucose, Hemin and Pen-Strep stock solutions)

Cat.-No.	Components	Size
ML-432S	20 g YS-1 powder 15 g YS-2 powder 5 ml Glucose 5 ml Pen-Strep 2 ml Hemin	For 1 L medium
ML-432L	100 g YS-1 powder 75 g YS-2 powder 25 ml Glucose 25 ml Pen-Strep 10 ml Hemin	For 5 L medium
ML-432XL	200 g YS-1 powder 150 g YS-2 powder 50 ml Glucose 50 ml Pen-Strep 20 ml Hemin	For 10 L medium
ML-432XXL	1000 g YS-1 powder 750 g YS-2 powder 250 ml Glucose 250 ml Pen-Strep 100 ml Hemin	For 50 L medium

Description:

Five-component Kit for preparation of complex *Leishmania* cultivation medium LEXSY YS.

The kit contains:

LEXSY YS-1 and YS-2, powder

200x Glucose stock solution:

20% solution of glucose in aqua dest. sterile filtrated

500x Hemin stock solution (ML-104):

0,25% solution of porcine Hemin in 50% Triethanolamine, sterile filtrated

200x Pen-Strep stock solution (ML-105):

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 filtrated

Storage conditions:

LEXSY YS-1 and YS-2 powder:

Store at room temperature
 Stable for 3 years

Glucose stock solution:

Store at 4°C in the dark
 Stable for 6 month

Hemin stock solution:

Store at 4°C in the dark
 Stable for 12 month

Pen-Strep stock solution:

Store at -20°C
 Stable for 12 month.

Preparation of complete *Leishmania* cultivation medium LEXSY YS:

Dissolve 20 g YS-1 powder and 15 g YS-2 powder in 500 ml aqua dest. each.

Autoclave YS-1 for 20 min and YS-2 for 15 min at 121°C in batches of 0.5 or 1 litre. Control temperature and time of sterilization with a reference with the same volume, since overexposure leads to decomposition of ingredients and a dark brown color of the medium, which adversely affects growth of LEXSY strains. Keep at room temperature. Stable for 6 month.

Prior to use mix YS-1 and YS-2 at 1:1 ratio and add to 1 L preformulated LEXSY YS:

5.0 ml of 200x Glucose solution (1 g/L)

5.0 ml of 200x Pen-Strep solution

2.0 ml of 500x Hemin stock solution to a final concentration of 5 µg Hemin /ml

Final pH = 7.2 ± 0.2

Store in the dark at 4°C

Completed medium is useable for up to 2 weeks.

Find more information in the User's Guide for LEXSY Expression Kits. Download pdf from the LEXSY section of our website

<http://www.jenabioscience.com/LEXSY>

FOR RESEARCH USE ONLY.

NOT INTENDED FOR HUMAN OR ANIMAL DIAGNOSTIC OR THERAPEUTIC USES.