

**LEXSY Neo**

sterile ready-to-go stock solution
Geneticin, G-418

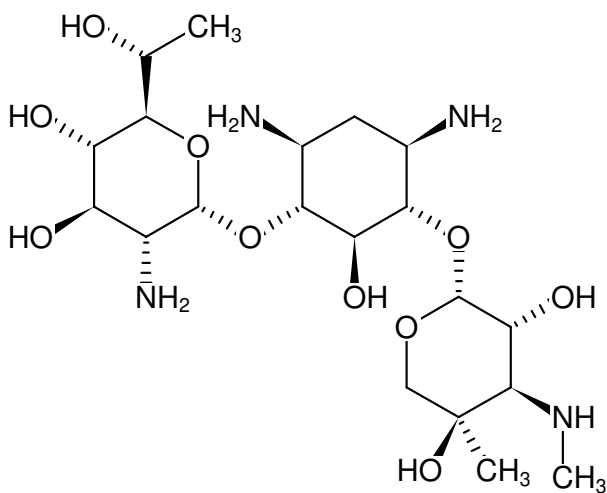
Cat. No.	Amount
AB-105S	1 ml (50 mg/ml)
AB-105L	5 x 1 ml (50 mg/ml)

Description:

LEXSY Neo is an aminoglycoside antibiotic used for the selection of recombinant *Leishmania* strains. It can also be used for the selection of other recombinant organisms expressing the APH gene encoding aminoglycoside 3'-phosphotransferase.

Selection:

For selection of recombinant *Leishmania* strains LEXSY Neo is added to the growth medium to a final concentration of 50 µg/ml.



Structural formula of LEXSY Neo

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

Shipping: shipped on gel packs

Storage Conditions: Store at -20 °C

Shelf Life: 12 months

Molecular Formula: C₂₀H₄₀N₄O₁₀

Molecular Weight: 496.6

Form: liquid

Color: colorless

Concentration: 50 mg/ml