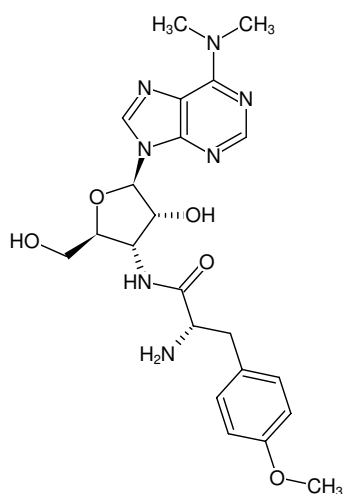


**LEXSY Puro**

sterile ready-to-go stock solution
Puromycin

Cat. No.	Amount
AB-108L	5 x 1 ml (10 mg/ml)



Structural formula of LEXSY Puro

For research use only! 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: C₂₂H₂₉N₇O₅

Molecular Weight: 471.5 g/mol

CAS#: 58-58-2

EC number: 200-387-8

Form: liquid

Color: colorless

Concentration: 10 mg/ml

Description:

LEXSY Puro is a nucleoside antibiotic used for the selection of recombinant *Leishmania* strains. It can also be used for the selection of other recombinant organisms expressing the *puro* marker gene (puromycin N-acetyltransferase from *Streptomyces alboniger*).

Selection:

For selection of recombinant *Leishmania* strains LEXSY Puro is added to the growth medium to a final concentration of 10 µg/ml.