

**LEXSY Bleo**

sterile ready-to-go stock solution
Zeocin, Phleomycin D1

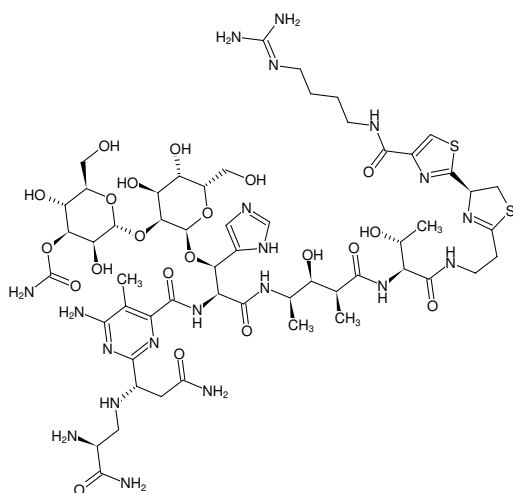
Cat. No.	Amount
AB-103S	1 ml (100 mg/ml)
AB-103L	5 ml (100 mg/ml)

Description:

LEXSY Bleo is a glycopeptide antibiotic used for the selection of recombinant *Leishmania* strains. It can also be used for the selection of other recombinant organisms expressing the *ble* gene encoding bleomycin resistance protein from *S. hindustans*.

Selection:

For selection of recombinant *Leishmania* strains LEXSY Bleo is added to growth medium to a final concentration of 100 µg/ml.



Structural formula of LEXSY Bleo

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: C₅₅H₈₆N₂₀O₂₁S₂

Molecular Weight: 1427.5 g/mol

Form: aqueous solution

Color: blue

Concentration: 100 mg/ml

pH: 7.25