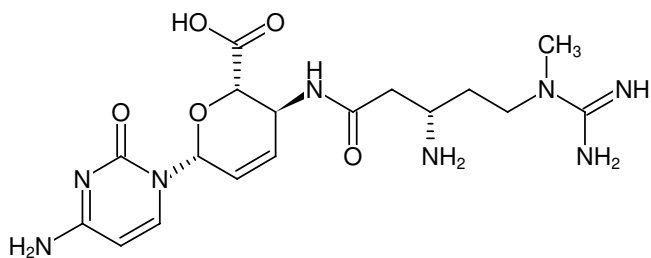


**LEXSY BSD**

sterile ready-to-go stock solution
Blasticidin S

Cat. No.	Amount
AB-107L	5 x 1 ml (10 mg/ml)



Structural formula of LEXSY BSD

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: C₁₇H₂₆N₈O₅

Molecular Weight: 422.4 g/mol

CAS#: 3513-03-9

Form: liquid

Color: colorless

Concentration: 10 mg/ml

Description:

Blasticidin 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 *bsd* marker gene (blasticidin-resistance of *Aspergillus terreus*).

Selection:

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