

**LEXSY Tet**

sterile ready-to-go stock solution
(for inducible protein expression in **LEXSY host T7-TR**)
Tetracycline

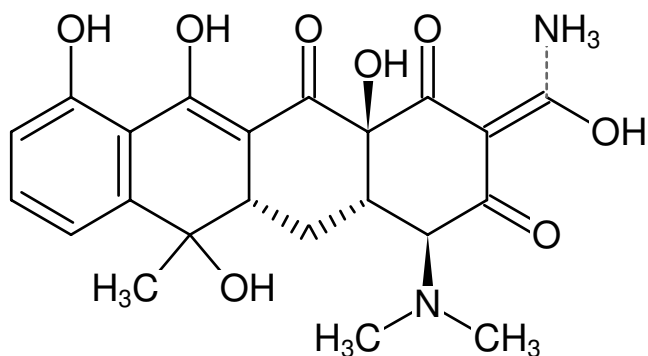
Cat. No.	Amount
AB-106S	5 ml (10 mg/ml)
AB-106L	10 ml (10 mg/ml)
AB-106XL	50 ml (10 mg/ml)

Description:

Tetracycline is an antibiotic which is used at sublethal concentrations for the induction of protein expression in the *Leishmania tarentolae* T7-TR strain.

Induction:

For induction of gene expression in recombinant *Leishmania* T7-TR strains LEXSY Tet is added to the growth medium to a final concentration of 10 µg/ml.



Structural formula of LEXSY Tet

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

Additional Storage Conditions: It is recommended to aliquot the content of the vial to avoid frequent freeze-thaw cycles. Turbidity of solution upon thawing does not compromise induction and is not a sign of malfunction. Mix well before use.

Shelf Life: 12 months

Molecular Formula: C₂₂H₂₄N₂O₈

Molecular Weight: 444.4

CAS#: 64-75-5

EC number: 200-593-8

Form: suspension

Color: yellow

Concentration: 10 mg/ml