

LEXSY-Tet (Tetracycline)

ready-to-use 1.000x stock solution (10 mg/ml)

	Cat.-No.	Size
	AB-106S	1 ml for 1 L
	AB-106L	5 ml for 5 L
	AB-106XL	10 ml for 10 L
	AB-106XXL	50 ml for 50 L

Description:

Contains 10 mg/ml Tetracycline antibiotic in water.
Sterile 1.000x stock solution.
For induction of gene expression in *Leishmania tarentolae* T7-TR strains.

Storage conditions:

Store at -20°C.
Stable for 12 month.
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.

Induction:

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

Please, find detailed information on induction conditions in the User's Guide for the LEXSY Expression Kit for inducible gene expression (download pdf from the LEXSY section of our website).

**FOR RESEARCH USE ONLY.
NOT INTENDED FOR HUMAN OR ANIMAL
DIAGNOSTIC OR THERAPEUTIC USES.**