

## LEXSY NTC

ready-to-use stock solution (100 mg/ml)

	Cat.-No.	Amount
	AB-101S	1 ml
	AB-101L	5 ml
	AB-101-10ML	10 ml
	AB-101-50ML	50 ml

### Field of use:

Streptothricin antibiotic Nourseothricin (NTC) for a broad spectrum of bacteria and other unicellular or complex organisms. Preferred selection antibiotic for genetically modified Gram-positive and Gram-negative bacteria, yeast, filamentous fungi, protozoa, microalgae, plants and many more. Selection of recombinant strains is based on inactivation of NTC by mono-acetylation of  $\beta$ -amino group of the  $\beta$ -lysine by Nourseothricin N-acetyltransferase the product of the *sat1* or *nat1* genes.

### Description:

Contains 100 mg/ml of sterile filtrated Nourseothricin in water.

### Storage conditions:

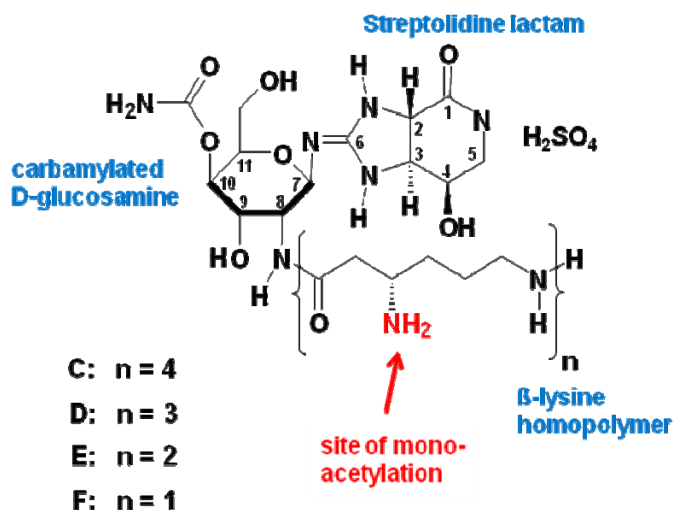
Store at -20°C.

Stable for 12 months.

### Selection:

For selection of recombinant *Leishmania* strains LEXSY NTC is added to growth medium to a final concentration of 100  $\mu$ g/ml.

**For selection of other species  
 please refer to Table 1**



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

## LEXSY NTC

### ready-to-use stock solution (100 mg/ml)

Table 1

Group	Species	MIC* µg/ml	Selection conc. µg/ml
Gram-negative bacteria	<i>Agrobacterium tumefaciens</i>		100
	<i>Escherichia coli</i>	2-12	50
	<i>Francisella tularensis</i>		50
	<i>Pseudomonas aeruginosa</i>	50	100
Gram-positive bacteria	<i>Bacillus subtilis</i>	5	50
	<i>Enterococcus faecium</i>	8-256	500
	<i>Staphylococcus aureus</i>	2-12	50
Streptomyces	<i>Streptomyces lividans</i>	6	100
Yeast	<i>Candida albicans</i>	200	250-450
	<i>Hansenula polymorpha</i>		100
	<i>Kluyveromyces lactis</i>		50
	<i>Pichia pastoris</i>		100
	<i>Saccharomyces cerevisiae</i>	25	75-100
	<i>Schizosaccharomyces pombe</i>	40	100
Other Ascomycota	<i>Acremonium chrysogenum</i>		25
	<i>Aspergillus nidulans</i>		120
	<i>Cryphonectria parasitica</i>		100
	<i>Neurospora crassa</i>		200
	<i>Penicillium chrysogenum</i>		150-200
	<i>Podospora anserina</i>		50
	<i>Sordaria macrospora</i>		50
	<i>Trichophyton mentagrophytes</i>		50
Basidiomycota	<i>Cryptococcus neoformans</i>		100
	<i>Schizophyllum commune</i>	3	8
	<i>Ustilago maydis</i>		75-100
Protozoa	<i>Leishmania tarentolae</i> , major etc.		100
	<i>Phytomonas serpens</i>		100
	<i>Plasmodium falciparum</i>	75** (IC50)	
	<i>Toxoplasma gondii</i>		500
Microalgae	<i>Phaeodactylum tricornutum</i>		50-250
	<i>Thalassiosira pseudonana</i>		100
Plants	<i>Arabidopsis thaliana</i>	20	50-200
	<i>Daucus carota</i>		100
	<i>Lotus corniculatus</i>		50
	<i>Nicotiana tabacum</i>		100
	<i>Oryza sativa</i>	20	200

## LEXSY NTC

### ready-to-use stock solution (100 mg/ml)

#### Selected References

- Alshahni *et al.* (2010) Nourseothricin acetyltransferase: a new dominant selectable marker for the dermatophyte *Trichophyton mentagrophytes*. *Medical Mycology* **48**:665
- Cox *et al.* (2003) Superoxide Dismutase Influences the Virulence of *Cryptococcus neoformans* by Affecting Growth within Macrophages. *Infect. Immunity* **71**: 173
- EI-Khoury *et al.* (2008) Gene deletion and allelic replacement in the filamentous fungus *Podospora anserina*. *Curr Genet* **53**: 249
- Gold *et al.* (1994) Three selectable markers for transformation of *Ustilago maydis*. *Gene* **142**: 225
- Goldstein *et al.* (1999) Three New Dominant Drug Resistance Cassettes for Gene Disruption in *Saccharomyces cerevisiae*. *Yeast* **15**: 1541
- Hamano *et al.* (2006) A Novel Enzyme Conferring Streptothricin Resistance Alters the Toxicity of Streptothricin D from Broad-spectrum to Bacteria-specific. *J. Biol. Chem.* **281**: 16842
- Hentges *et al.* (2005) Three novel antibiotic marker cassettes for gene disruption and marker switching in *Schizosaccharomyces pombe*. *Yeast* **22**: 1013
- Hoff *et al.* (2009) Homologous recombination in the antibiotic producer *Penicillium chrysogenum*: strain ΔPcku70 shows up-regulation of genes from the HOG pathway. *Appl Microbiol Biotechnol* **85**:1081
- Jacobs *et al.* (2009) Engineering complex-type N-glycosylation in *Pichia pastoris* using GlycoSwitch technology. *Nature Protocols* **4**: 58
- Jelenska *et al.* (2000) Streptothricin resistance as a novel selectable marker for transgenic plant cells. *Plant Cell Reports* **19**: 298
- Ji *et al.* (2007) Two New Members of Streptothricin Class Antibiotics from *Streptomyces qinlingensis* sp. Nov. *J. Antibiot.* **60**: 739
- Joshi *et al.* (1995) The gene encoding streptothricin acetyltransferase (*sat*) as a selectable marker for *Leishmania* expression vectors. *Gene* **156**: 145
- Kloti *et al.* (2004) Selectable marker in plants. *United States Patent* 6696621
- Kojic *et al.* (2000) Shuttle vectors for genetic manipulations in *Ustilago maydis*. *Can. J. Microbiology* **46**: 333
- Kooistra *et al.* (2004) Efficient gene targeting in *Kluyveromyces lactis*. *Yeast* **21**: 781
- Kück *et al.* (2006) Application of the nourseothricin acetyltransferase gene (*nat1*) as dominant marker for the transformation of filamentous fungi. *Fungal Genetics Newsletter* **53**: 9
- Laubinger (2008) Analysis of Neuronal Diseases in the Model Organism *Aspergillus nidulans*. *Dissertation Univ. Göttingen*
- Lukeš *et al.* (2006) Translational initiation in *Leishmania tarentolae* and *Phytomonas serpens* (Kinetoplastida) is strongly influenced by pre-ATG triplet and its 5' sequence context. *Molecular & Biochemical Parasitology* **148**: 125
- Maier *et al.* (2006) In Vivo *Himar1*-Based Transposon Mutagenesis of *Francisella tularensis*. *Appl. Environment. Microbiol.* **72**: 1878
- McDade *et al.* (2001) A new dominant selectable marker for use in *Cryptococcus neoformans*. *Med. Mycol.* **39**: 151
- Mamoun *et al.* (1999) A set of independent selectable markers for transfection of the human malaria parasite *Plasmodium falciparum*. *PNAS* **96**: 8716
- Mandal *et al.* (2009) High-throughput screening of amastigotes of *Leishmania donovani* clinical isolates against drugs using colorimetric β-lactamase assay. *Ind. J. Exptl. Biol.* **47**: 475
- Nagotu *et al.* (2008) Peroxisome Fission in *Hansenula polymorpha* Requires Mdv1 and Fis1, Two Proteins Also Involved in Mitochondrial Fission. *Traffic* **9**: 1471
- Ohm *et al.* (2010) An efficient gene deletion procedure for the mushroom-forming basidiomycete *Schizophyllum commune*. *World J Microbiol Biotechnol* **26**:1919
- Poulsen *et al.* (2006) Molecular genetic manipulation of the diatom *Thalassiosira pseudonana* (bacillariophyceae) *Phycology* **42**: 1059
- Shen *et al.* (2005) CaNAT1, a heterologous dominant selectable marker for transformation of *Candida albicans* and other pathogenic *Candida* species. *Infection and Immunity*, **73**: 1239
- Smith *et al.* (2007) Two yeast plasmids that confer nourseothricin-dihydrogen sulfate and hygromycin B resistance in *Neurospora crassa* and *Cryphonectria parasitica*. *Fungal Genetics Newsletter* **54**: 12
- Van *et al.* (2006) Nourseothricin Acetyltransferase: A Positive Selectable Marker for *Toxoplasma gondii*. *J. Parasitol.* **92**: 668
- van Peer *et al.* (2009) Phleomycin Increases transformation efficiency and promotes single integrations in *Schizophyllum commune*. *Appl. Environ. Microbiol.* **75**, 1243-1247
- Werner *et al.* (2001) Aminoglycoside-Streptothricin Resistance Gene Cluster *aadE-sat4-aphA-3* Disseminated among Multiresistant Isolates of *Enterococcus faecium*. *Antimicrob. Agents Chemotherapy* **45**: 3267
- Zaslavskaja *et al.* 2000 Transformation of the diatom *Phaeodactylum tricoratum* (bacillariophyceae) with a variety of selectable marker and reporter genes. *J. Phycol.* **36**: 379

\* MIC: Minimal inhibitory concentration

\*\* IC50: Concentration inhibiting growth by 50%