



# Nourseothricin past, present and future

Reinhard Breitling

Jena Bioscience GmbH  
Löbstedter Str. 80  
07749 Jena, Germany  
Tel.: +49-3641-628-5000  
Fax: +49-3641-628-5100

e-Mail: [info@jenabioscience.com](mailto:info@jenabioscience.com)

# Jena Bioscience provides >5000 life science products

**Jena Bioscience**

Products & Ordering | Order Information | About us | Contact | Downloads | Blog

Search

IFTA AG  
Certified QMS according to  
DIN EN ISO 9001  
Reg. No. IC 03214 034

**News**

May 17, 2011 **New in Click Chemistry: Nucleotide Azides**  
We have completed our [Click Chemistry catalog section](#) by a selection of [Nucleotide Azides](#) for labeling and detection using "Click" reactions.

Apr 28, 2011 **Pi sampling strategy**  
The new [Pi-minimal Screen](#) was developed at the MRC Cambridge. The incomplete factorial design assures diversity amongst the crystallization conditions ideal for initial screening.

Apr 28, 2011 **Pi sampling applied to membrane proteins**  
Use our new [Pi-PEG Screen](#) and benefit from the experience of membrane protein crystallization at the MRC Cambridge!

Apr 21, 2011 **Maximum durability and rigidity combined with lowest X-ray background!**  
The new [Dual-Thickness MicroMounts™](#) are now available!

Apr 20, 2011 **New Click Chemistry Reagents**  
We have again extended our [Click Chemistry section](#) by [Alkyne-containing Phosphoramidites](#) and [CPGs](#) as well as by several new [Azides of Fluorescent Dyes!](#)

**Nucleosides, Nucleotides and their Analogs**

**Molecular Biology**

**Fluorescent Probes**

**Macromolecular Crystallography**

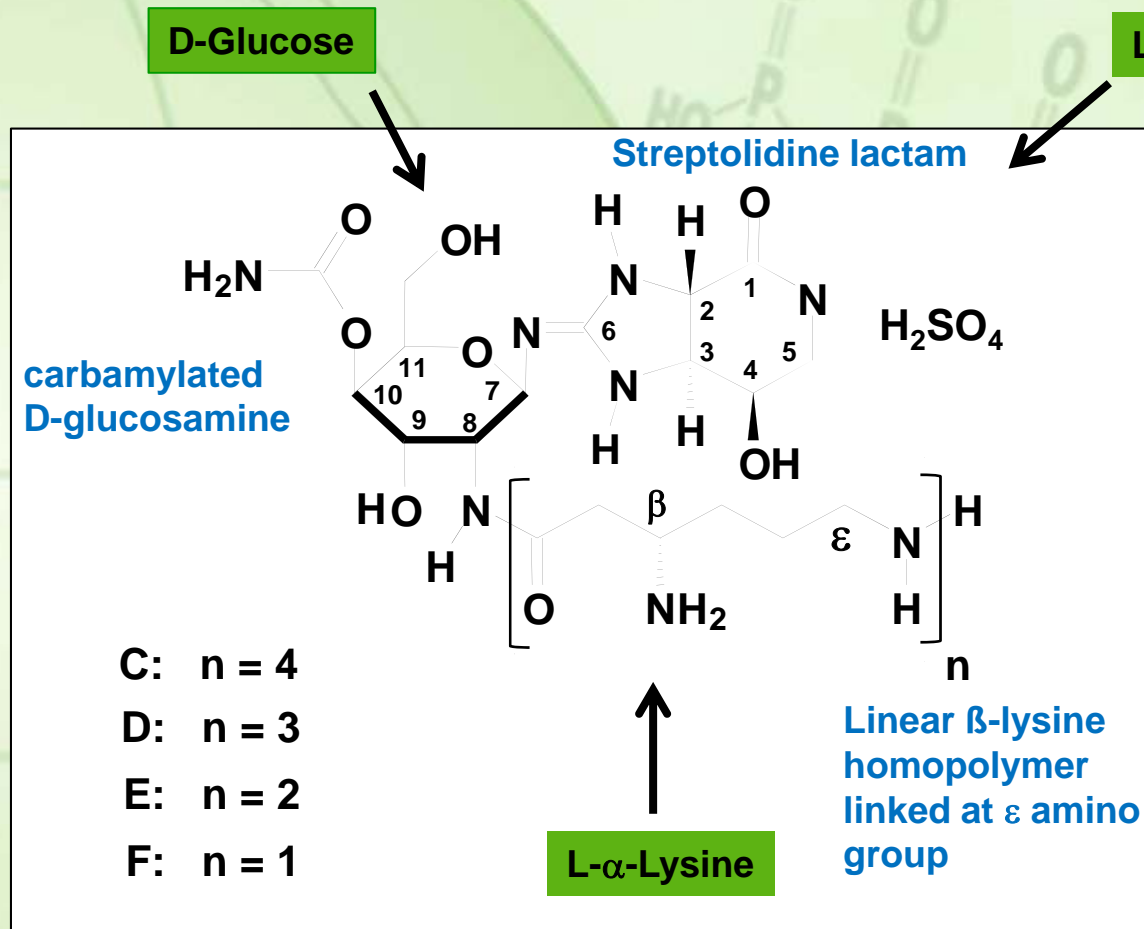
**Eukaryotic Expression System LEXSY**

**Recombinant Proteins**

**Biochemistry**

**Affinity Chromatography**

# Nourseothricin is a aminoglycoside glycopeptide (nucleoside peptide) antibiotic of the Streptothricin class



Secondary metabolite of Nystatin producer *Streptomyces noursei* ZIMET JA3890 (ATCC 11455) (Bradler *et al.* 1963)

Synthesised from Glucose, Arginine and Lysine by convergent pathways (Jackson *et al.* 2002 and ref.)

Natural mixture of streptothricins C, D, E & F; D + F >85%

Substance produced exclusively at HKI Jena



## Members of the Streptothricin class antibiotics differ by the number of $\beta$ -lysine residues (n = 1-7)

Producer	Antibiotic	Reference	
<i>Streptomyces lavendulae</i>	Streptothricin F	Waksman <i>et al.</i> 1942	F n=1
<i>Streptomyces griseus</i>	Streptothricin C, D, E, F (C + F > D + E)	Reynolds <i>et al.</i> 1947	E n=2
<i>Streptomyces noursei</i>	Streptothricin C, D, E, F (D + F >85%)	Bradler <i>et al.</i> 1963	D n=3
<i>Streptomyces rochei</i>	Streptothricin F , D, E	Singh <i>et al.</i> 1983	C n=4
<i>Streptomyces spp.</i> SNUS 8810-111	Streptothricin D + mod	Kim <i>et al.</i> 1994	B n=5
<i>Streptomyces quinlingensis</i> <i>sp.nov.</i>	Streptothricin D, F + mod	Ji <i>et al.</i> 2007	A n=6
			X n=7

- First member: Streptothricin F (n=1)
- Proposal of structural formula (van Tamelen *et al.* 1961)
- Separation of Streptothricins by ion exchange chromatography (Reshetov *et al.* 1964)
- Chemical structure and general formula (n= 1-7) (Khokhlov *et al.* 1964-1978)
- Total chemical structure (Kusumoto *et al.* 1982)

→ Streptolin, Geomycin, Phytobacteriomycin, Racemomycin, Pleocidin, Polymycin, Yazumycin, Grisein, Nourseothricin ... are mixtures of Streptothricins

# Nourseothricin (NTC) was applied as an ergotropic agent

- Broad spectrum antibacterial effect
- Not used for therapeutic purposes (human or veterinary) because of nephrotoxicity
- Not resorbed by intestinum wall
- More efficient biomass production in animal farms due to inhibition of growth of (competing) microflora of digestive tract



- Produced at large scale as bentonite adsorbate of mycelium by Jenapharm in Jena
- Controlled application at selected sites in Germany in the 1980th



## **Streptothricin resistances were found at sites where Nourseothricin was feeded to pigs**

- **Resistance plasmids were found in *E. coli* from pigs and from employees in pig farms and their family members (Hummel *et al.* 1986)**
- **Resistance plasmids were also found in man without contact to animal farms but living in territories where Nourseothricin was applied as ergotropic agent (Hummel *et al.* 1986)**
- **The resistance plasmids found in *E. coli* from man were similar to the plasmids of *E. coli* from pigs and were of different incompatibility groups (Hummel *et al.* 1986)**
- **Hybridization to bacterial Streptothricin resistance gene probes was also observed with plasmids isolated a long time before the application of streptothricins (Tietze *et al.* 1990)**
- **The streptothricin resistance determinants were found to be linked to other resistance genes like streptomycin/spectinomycin- and trimethoprim-resistances on bacterial transposons (Sundström *et al.* 1991).**

## Three bacterial Streptothricin resistance genes where found at sites related to ergotropic use of streptothricins

Gene	Source	Reference	Linkage
sat1*	Bacterial transposon Tn 1825	Heim <i>et al.</i> 1989	<i>sat - aadA1</i>
sat2*	Bacterial transposon Tn 1826	Tietze <i>et al.</i> 1990a	<i>sat - aadA1</i>
sat*	Bacterial transposon Tn7	Sundström <i>et al.</i> 1991	<i>dhfrAI - sat - aadA1</i>
sat3	E. coli Plasmids pIE636, pIE637 and pIE639	Seltmann 1985 Tietze <i>et al.</i> 1990b	
sat4	<i>Campylobacter coli</i> BE/G4 <i>Enterococcus faecium</i> <i>S. aureus</i> Tn 5405	Jacob <i>et al.</i> 1994 Werner <i>et al.</i> 2001 Debrise <i>et al.</i> 1996	<i>aadE - sat4 - aphA-3</i>

sat1 = sat2 = sat Tn7 (Sundström *et al.* 1991)

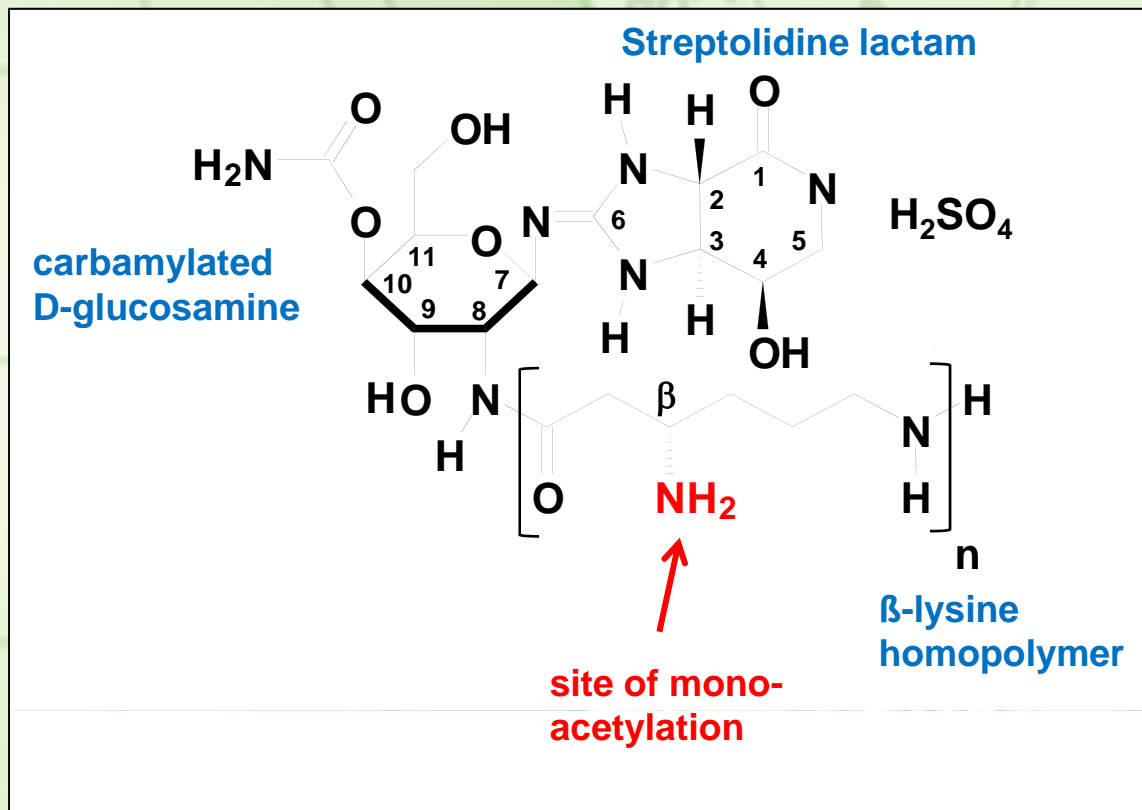
*aadA1* = streptomycin/spectinomycin resistance

*aadE* = aminoglycoside resistance

*aphA-3* = kanamycin resistance

*dhfrAI* = trimethoprim resistance

## The *sat* resistance genes encode a streptothricin N-acetyltransferase



Streptothricins are inactivated by monoacetylation of the β-amino group of the β-lysine residue

The 524 bp *sat* gene encodes a small protein of 174 aa (20 kDa)

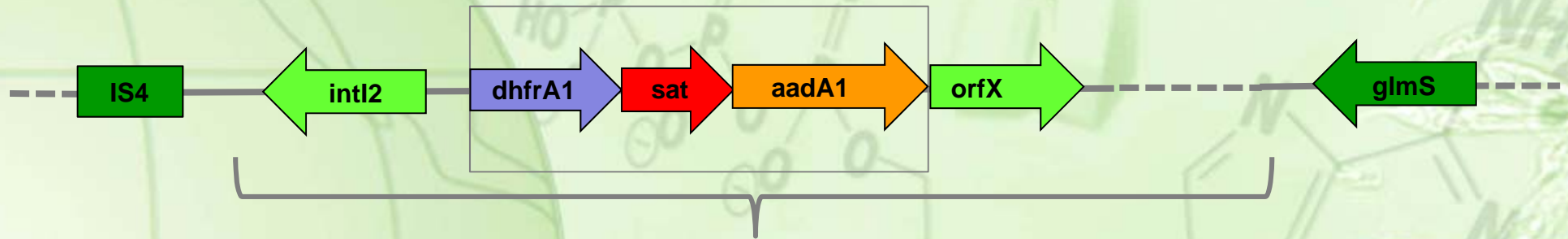


## Sat genes were found worldwide also in numerous other bacteria

Species	Source	Reference	Linkage
<i>Aerococcus viridans</i>	animal site	Byrne-Bailey <i>et al.</i> 2008	Class1/2 integron
<i>Acinetobacter baumannii</i>		Ploy <i>et al.</i> 2000, Ramirez <i>et al.</i> 2005	Class 2 integron Tn7::In2-8
<i>Burkholderia cenocepacia</i>	clinical isolate	Ramirez <i>et al.</i> 2005	Class 2 integron Tn7::In2-1
<i>Citrobacter freundii</i>	beef cuttle AU	Barlow <i>et al.</i> 2006	Class 2 integron
<i>Enterobacter cloacae</i>	clinical isolate	Ramirez <i>et al.</i> 2005	Class2 integron with IS1 (Tn7::In2-10)
<i>Enterococcus faecalis</i>	clinical isolate	Xu <i>et al.</i> 2010	Class2 integron
<i>Klebsiella oxytoca &amp; pneumonia</i>	clinical isolate	Xu <i>et al.</i> 2007	Class2 integron
<i>Morganella morganii</i>	clinical isolate	Power <i>et al.</i> 2005	Class 2 integron
<i>Proteus mirabilis</i>	meat products	Kim <i>et al.</i> 2004	
<i>Pseudomonas aeruginosa</i> Tn7		Ramirez <i>et al.</i> 2005	MDR
<i>Psychrobacter maritimus &amp; sp.</i>	animal site	Byrne-Bailey <i>et al.</i> 2008	Class1/2 integron
<i>Raoultella terrigena</i>	clinical isolate	Ramirez <i>et al.</i> 2005	Class2 integron-Tn7
<i>Salmonella enterica</i> SV enteritidis	clinical isolate	Ahmed <i>et al.</i> 2005	Class 2 integron
<i>Serratia marcescens</i>	clinical isolate	Crowley <i>et al.</i> 2006	Class2 integron
<i>Shigella flexneri &amp; sonnei</i>	clinical isolate	Halloran <i>et al.</i> 2002, Pan <i>et al.</i> 2006	Class1/2 integron MDR
<i>Vibrio cholerae</i>	clinical isolate	Coelho <i>et al.</i> 1995, Ahmed <i>et al.</i> 2006	Class1/2 integron

Protein Blast "sat": 99-100% identity

## The integrons contained multiple antibiotic resistance markers



Tn7 containing a class 2 integron (14 kbp) isolated from *Shigella sonnei* after Pan *et al.* 2006

- Co-selection of resistances to therapeutic antibiotics (Streptomycin)
- Ergotropic use of Nourseothricin was stopped after 1989 in Germany

## Streptothricin producers harbour self-resistance genes encoding a streptothricin N-acetyltransferase

Gene	Species	Reference
<i>stat</i>	<i>Streptomyces lavendulae</i>	Horinouchi et al. 1987
<i>nat1 &amp; 2</i>	<i>Streptomyces noursei</i>	Krügel et al. 1993
<i>gsr</i>	<i>Streptomyces griseus</i>	Sezonov et al. 1990
<i>sttR</i>	<i>Streptomyces rochei</i>	Anukool et al. 2004
<i>NAT_SF</i>	<i>Streptomyces ambofaciens</i>	Choulet et al. 2006
<i>NAT_SF</i>	<i>Streptomyces pactum</i>	Ito et al. 2008
<i>NAT_SF</i>	<i>Streptomyces sp. C</i>	Fischbach et al. 2009
<i>NAT_SF</i>	<i>Streptomyces sp. e14</i>	Fischbach et al. 2010
<i>NAT_SF</i>	<i>Streptomyces roseosporus</i>	genome shotgun sequences 2010
<i>NAT_SF</i>	<i>Streptomyces cattleya</i>	Centre National de Sequencage 2011

Protein Blast “*nat1*”: 66-78% similarity

→ opens the way for genetic manipulation of sensitive organisms in combination with streptothricin antibiotics



# Nourseothricin: A superior selection antibiotic in molecular genetics

## Field of use

- Extraordinarily broad spectrum of sensitive bacteria and eukaryotic organisms
- Excellent selection antibiotic for genetic modification of
  - Gram-positive and Gram-negative bacteria
  - Yeast and filamentous fungi
  - Protozoa and microalgae
  - Plants ... and more

## Mechanism of Action

- Antibiotic effect of Nourseothricin through inhibition of protein biosynthesis and induction of miscoding
- Resistance to Nourseothricin conferred by *sat* or *nat* marker genes
- Product of the resistance gene - Nourseothricin N-acetyltransferase - inactivates NTC by monoacetylation of  $\beta$ -amino group of the  $\beta$ -lysine residue

## Advantages

- Low or no background: Resistance protein is localized intracellularly and cannot be degraded in the cell culture medium
- Not used in human or veterinary medicine, therefore, no conflict with regulatory requirements
- No cross-reactivity with other aminoglycosid antibiotics such as Hygromycin or Geneticin
- No cross-resistance with therapeutic antibiotics
- Long-term stable as powder or solution
- Highly soluble in water (1 g/L)

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, major etc.</i>	30-50	100
	<i>Phytomonas serpens</i>		100
	<i>Plasmodium falciparum</i>	75**	
	<i>Toxoplasma gondii</i>		500
Microalgae	<i>Phaeodactylum tricoratum</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

**Nourseothricin is used in about 50 recombinant host-vector systems**

**permanently increasing number of species for genetic engineering will further extend its application**

\*MIC:  
Minimal inhibitory concentration

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

# Acknowledgments

**Dr. Walter Werner & Gisela Werner  
WERNER BioAgents, Jena**

**Dr. Hans Krügel  
Leibniz Institute Nat. Prod. Res. & Infect. Biology (HKI)**

**Thank you for your attention!**