

Staphylococcus aureus, *S. aureus* 8325-4 (Xen8.1)

Product Number: 119239

Material Provided: 1 Agar Plate

Storage Conditions: -80°C

Genetic Characteristics

Staphylococcus aureus-Xen8.1 was derived from the parental strain *S. aureus* 8325-4. *S. aureus*-Xen8.1 was engineered through transposition of Tn4001 luxABCDE on plasmid pXen-5. Xen8.1 possesses a single stable copy of the modified *P. luminescens* lux operon that was inserted in the δ -toxin coding region in the RNAIII transcript downstream of the agr P3 promoter on the bacterial chromosome.

Growth Characteristics

S. aureus-Xen8.1 grows well in Luria Bertani (LB) medium at 37°C under ambient aeration. *S. aureus*-Xen8.1 may also be grown selectively on LB or BHI agar containing 200µg/mL kanamycin.

Colonial Morphology

On LB agar, *S. aureus*-Xen8.1 appears as small (~1.5mm), cream-colored, opaque, smooth, circular colonies.

Growth Curve

S. aureus-Xen8.1 displays peak bioluminescence during early log-phase growth. Log-phase growth can be achieved after 1 to 1.5 hours of subculture in LB broth at 37°C, shaking at 150-200 rpm. For these broth culture conditions, an absorbance measurement at 600nm (against a LB blank) of 0.6 is roughly equivalent to 5×10^8 cfu/mL of *S. aureus*-Xen8.1 and the relative light intensity is 0.9 photons/sec/cell.

Virulence Factors

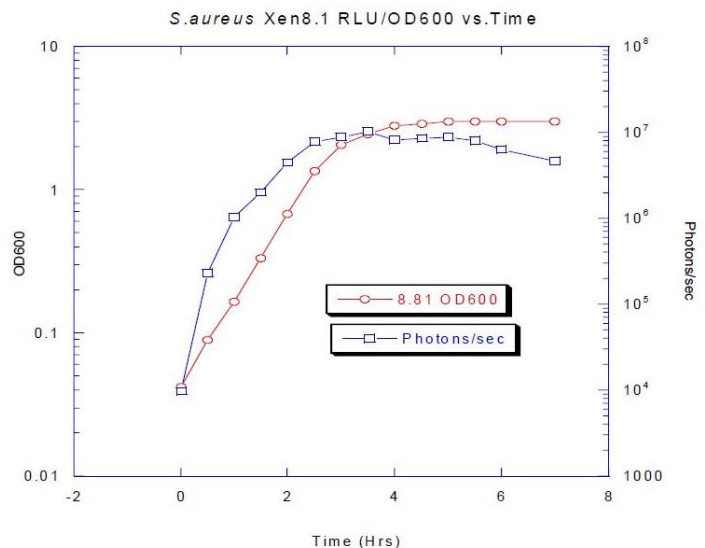
Hemolysins: β -hemolytic on TSA + 5% sheep blood

Capsule: literature cites that parental lacks a capsule (CP5 negative), cap5 mutant FEMS Microbiol Lett 1999 Jan 1;170(1):97-103.

DNase: Negative

NaCl: Tolerant via growth Mannitol Salts Agar

Coagulase: Positive in 24hrs



Biochemical Profile

A biochemical profile was obtained for *S. aureus*-Xen8.1 using the api 20 STAPH system available from bioMérieux.

Sugar Utilization	
D-Glucose	+
D-Fructose	+
D-Mannose	+
Maltose	+
Lactose	+
Trehalose	+
D-Mannitol	+
Xylitol	-
Raffinose	-
Xylose	-
D-Melibiose	+
Sucrose	+

Other Tests	
Nitrate Reduction	+
Alkaline Phosphatase	+
Voges Proskauer	-
□-methyl-D-glucoside	-
□-acetyl-glucoside	+
Arginine dihydrolase	+
Urease	+

Antibiotic Susceptibility

Disk Diffusion Data Disk diffusion tests were performed according to methods outlined in the NCCLS Approved Standard M2-A7.

Kirby-Bauer Disk Diffusion Test	
Sensitive to:	Resistant to:
Carbenicillin 100	Kanamycin 30
Gentamicin 20	
Penicillin G 10U	
Vancomycin 30	

MIC and MBC Data

MIC and MBC were determined using the macrodilution methods specified in the NCCLS Approved Standard M7-A5.

NCCLS Macrodilution MIC/MBC		
Antibiotic	MIC (□g/mL)	MBC (□g/mL)
Ceftriaxone	>32	NA
Ciprofloxacin	>8.0	NA
Erythromycin	1.0	1.0
Gentamicin	2.0-4.0	4.0-8.0
Kanamycin	>125	NA
Penicillin G	0.013-0.063	0.063-0.125
Tetracycline	0.25	0.5 (trailing endpoint)

The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

119239-R REV01

revvity