

Staphylococcus aureus, *S. aureus* UAMS-1 (Xen40)

Product Number: 119244

Material Provided: 1 Agar Plate Storage Conditions: -80°C

Genetic Characteristics

Staphylococcus aureus-Xen40 was derived from the parental strain *S. aureus* UAMS-1, a clinical isolate. *S. aureus*-Xen40 possesses a stable copy of the *Photorhabdus luminescens lux* operon at a single integration site on the bacterial chromosome.

Growth Characteristics

S. aureus-Xen40 grows well in Brain-Heart Infusion (BHI) broth at 37°C under ambient aeration. *S. aureus*-Xen40 may also be grown selectively on LB agar containing 200 µg/mL kanamycin.

Colonial Morphology

On LB agar, *S. aureus*-Xen40 appears as small (~1.5mm), cream-colored, opaque, smooth, circular colonies.

Growth Curve

Log-phase growth can be achieved after 1.5 to 3 hours of subculture in BHI broth at 37°C, shaking at 200 rpm. For the above broth culture conditions, an absorbance measurement at 600nm (against a LB blank) of 1 is roughly equivalent to 3x10⁸ cfu/mL of *S. aureus*-Xen40 and the relative light intensity is 0.013 photons/sec/cell.

Virulence Factors

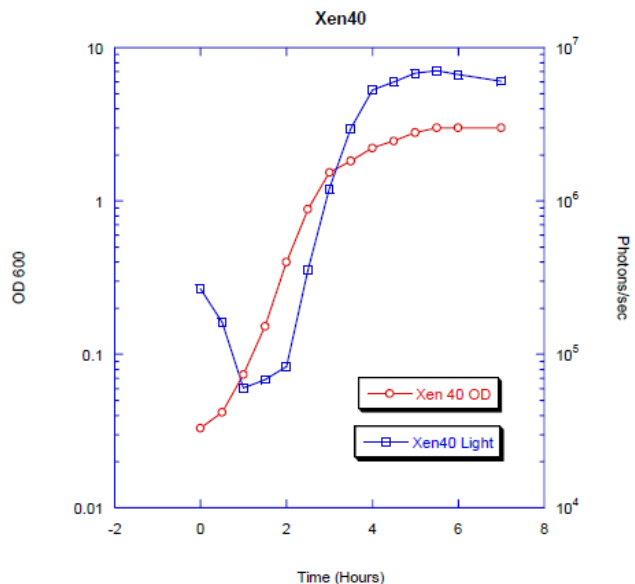
Hemolysins: β-hemolytic on TSA + 5% sheep blood

Coagulase: Positive in 24hrs (parental).

DNase: Positive

NaCl: Tolerant via growth Mannitol Salts Agar

Coagulase: Positive in 24hrs



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

119244-R REV01