

Shigella dysenteriae, S. dysenteriae 88A 6205 clinical isolate (Xen27)

Product Number: 119231

Material Provided: 1 Agar Plate

Storage Conditions: -80°C

Genetic Characteristics

Shigella dysenteriae-Xen27 was derived from the parental strain *S. dysenteriae* 88A 6205, a clinical isolate from California Department of Public Health. *S. dysenteriae*-Xen27 possesses a stable copy of the *Photorhabdus luminescens lux* operon on the bacterial chromosome.

Growth Characteristics

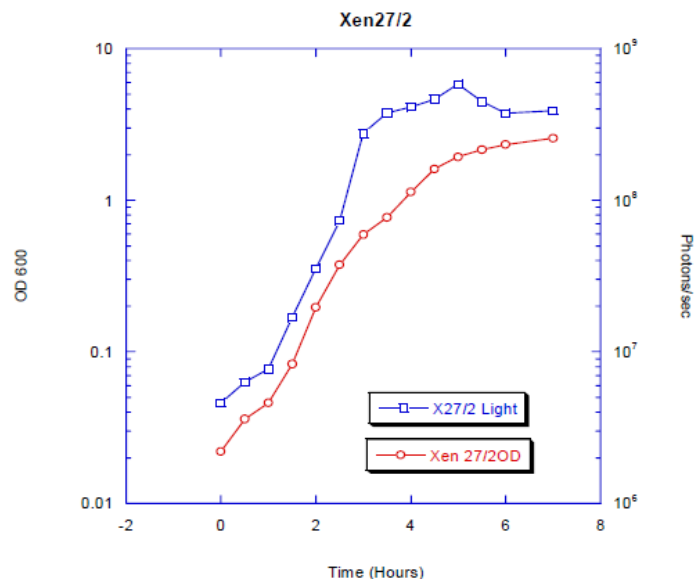
S. dysenteriae-Xen27 grows well in Luria Bertani (LB) medium at 37°C under ambient aeration. *S. dysenteriae*-Xen27 may also be grown selectively on LB agar containing 30µg/mL kanamycin.

Colonial Morphology

On LB plates, *S. dysenteriae*-Xen27 appears as small (1.5mm), cream, smooth, circular colonies after 24 hours incubation at 37°C.

Growth Curve

Log-phase growth can be achieved after 1 to 2.5 hours of subculture in LB broth at 37°C, shaking at 200 rpm. For the above broth culture conditions, an absorbance measurement at 600nm (against a LB blank) of 1.0 is roughly equivalent to 3.2×10^8 cfu/mL of *S. dysenteriae*-Xen27 and the relative light intensity is 10.3 photons/sec/cell.



Biochemical Profile

A biochemical profile was obtained for *P. aeruginosa* Xen 5 using the api 20NE system available from bioMérieux.

| Sugar Fermentation /Oxidation | | Other Tests | |
|-------------------------------|---|-----------------------------|---|
| Glucose | + | α-galactosidase | - |
| Mannitol | - | Arginine Dihydrolase | - |
| Inositol | - | Lysine Decarboxylase | - |
| Sorbitol | - | Ornithine Decarboxylase | - |
| Rhamnose | - | Citrate Utilization | - |
| Sucrose | - | H ₂ S Production | - |
| Melibiose | - | Urease | - |
| Amygdalin | - | Tryptophan Deaminase | - |
| Arabinose | - | Indole Production | - |
| | | Voges Proskauer | - |
| | | Gelatinase | - |
| | | Oxidase | - |

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: |
| Ampicillin 10 | |
| Carbenicillin 100 | |
| Chloramphenicol 30 | |
| Sulfamethoxazole/ Trimethoprim | |
| Tetracycline 30 | |

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

119231-R REV01