

Yersinia enterocolitica, Y. enterocolitica WS2589 (Xen25)

Product Number: 119233

Material Provided: 1 Agar Plate

Storage Conditions: -80°C

Genetic Characteristics

Yersinia enterocolitica-Xen25 was derived from the parental strain *Y. enterocolitica* WS2589, a clinical isolate from Weihenstephan Culture Collection. *Y. enterocolitica*-Xen25 possesses a stable copy of the *Photorhabdus luminescens lux* operon on the bacterial chromosome.

Growth Characteristics

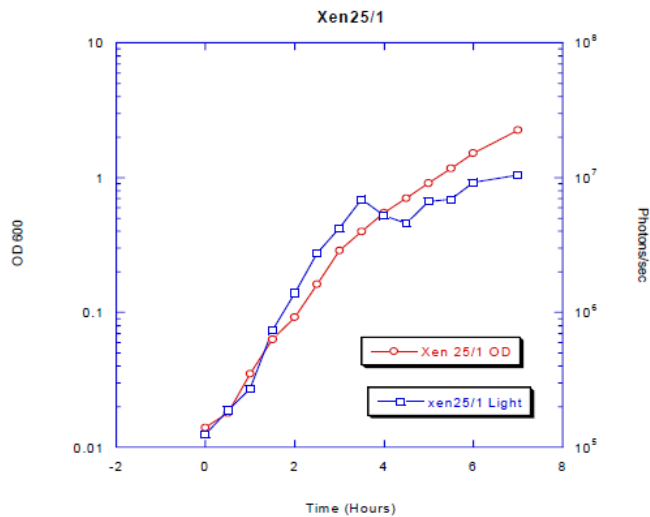
Y. enterocolitica-Xen25 grows well in Luria Bertani (LB) medium at 37°C under ambient aeration. *Y. enterocolitica*-Xen25 may also be grown selectively on LB agar containing 30µg/mL kanamycin.

Colonial Morphology

On LB plates, *Y. enterocolitica*-Xen25 appears as medium (2mm), cream colored, smooth, circular colonies after 24 hours incubation at 37°C.

Growth Curve

Log-phase growth can be achieved after 4 to 5 hours of subculture in BHI broth at 37°C, 5%CO₂, with no aeration in culture tubes. For the above broth culture conditions, an absorbance measurement at 600nm (against a BHI blank) of 0.5 is roughly equivalent to 2.4x10⁸ cfu/mL of *S. pneumoniae*-Xen11.



Biochemical Profile

A biochemical profile was obtained for *Yersinia enterocolitica*-Xen25 using the api 20E 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:
Chloramphenicol 30	Ampicillin 10
Sulfamethoxazole/ Trimethoprim	Carbenicillin 100
Tetracycline 30	Cephalothin 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

119233-R REV01

