

Listeria monocytogenes, *L. monocytogenes* ATCC 23074 (Xen19)

Product Number: 119237

Material Provided: 1 Agar Plate

Storage Conditions: -80°C

Genetic Characteristics

Listeria monocytogenes Xen 19 was derived from the parental strain *L. monocytogenes* ATCC23074, which was deposited by Stanford University as AT-25. *L.monocytogenes* Xen 19 possesses a stable copy of the modified *Photorhabdus luminescens lux* operon at a single integration site on the bacterial chromosome.

Growth Characteristics

L. monocytogenes Xen 19 grows well in Brain-Heart Infusion (BHI) broth at 37°C. It may also be grown on Trypticase Soy Agar (TSA). *L. monocytogenes* Xen 19 may be grown selectively on LB agar containing 200 mg/mL kanamycin.

Colonial Morphology

On TSA + 5% sheep blood plates, *L. monocytogenes* Xen 19 appears as small (approx 0.5mm), circular, gray colonies with large zones of beta hemolysis.

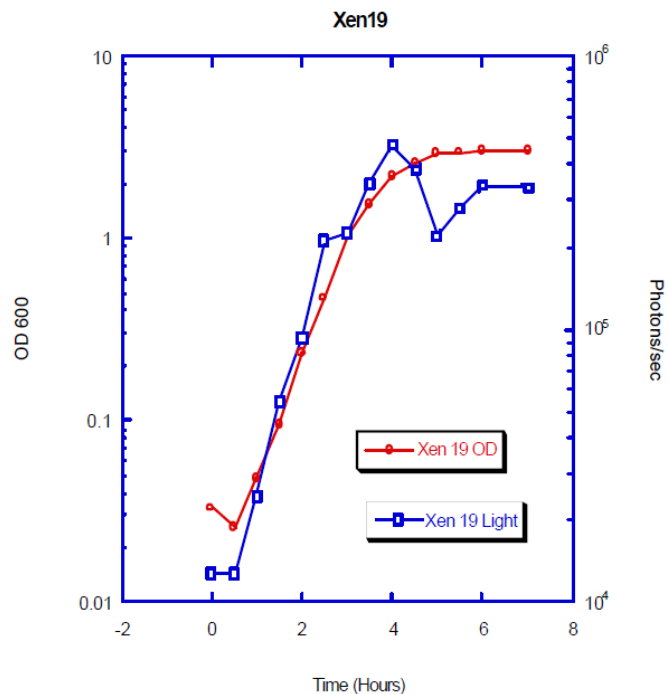
Growth Curve

Log-phase growth can be achieved after 1-3 hours of subculture in BHI broth at 37°C, shaking at 200 rpm. For these broth culture conditions, an absorbance measurement at 600 nm (against a BHI blank) of 1 is roughly equivalent to 4×10^8 cfu/ml of *L. monocytogenes* Xen 19 and the relative light intensity is 0.005 photons/sec/cell.

Virulence Factors

Hemolysins: \square -hemolytic on TSA + 5% sheep blood

Bacitracin: Susceptible



Biochemical Profile

A biochemical profile was obtained for *Listeria monocytogenes*-Xen19 using the api 20 Strep system available from bioMérieux.

Sugar Utilization	
Ribose	
Arabinose	
Mannitol	
Sorbitol	
Lactose	
Trehalose	
Inuline	
Raffinose	
Starch	
Glycogen	

Other Tests	
Voges Proskauer	
Hippurate	
Esculin	
Pyrrolidone Arylamidase	
<input type="checkbox"/> galactosidase	
<input type="checkbox"/> glucuronidase	
<input type="checkbox"/> galactosidase	
Alkaline Phosphatase	
Leucine Arylamidase	
Arginine Dihydrolase	

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

119237-R REV01

