

## Escherichia coli, E. coli WS2572 (Xen14)

Product Number: 119223

Material Provided: 1 Agar Plate

Storage Conditions: -80°C

### Genetic Characteristics

*Escherichia coli*-Xen14 was derived from the parental strain *Escherichia coli* WS2572, a clinical isolate from Weihenstephan Culture Collection. *E. coli*-Xen14 possesses a stable copy of the *Photorhabdus luminescens lux* operon on the bacterial chromosome.

### Growth Characteristics

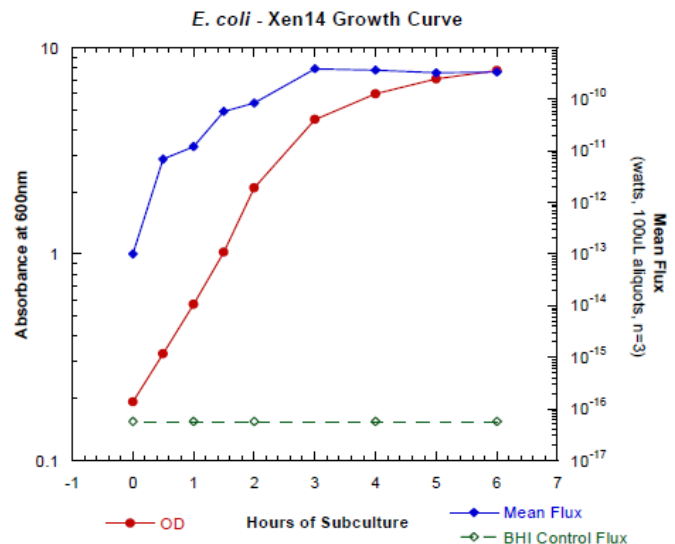
*E. coli*-Xen14 grows well in Luria Bertani (LB) medium at 37°C under ambient aeration. *E. coli*-Xen14 may also be grown selectively on LB agar containing 30µg/mL Kanamycin.

### Colonial Morphology

On LB plates, the *E. coli*-Xen14 appears as medium (~2mm), translucent, conical colonies after 24 hours incubation at 37°C.

### Growth Curve

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 the above broth culture conditions, an absorbance measurement at 600nm (against a LB blank) of 1.0 is roughly equivalent to  $1.1 \times 10^9$  cfu/mL of *Escherichia coli*-Xen14 and the relative light intensity is 1.01 photons/sec/cell.



## Biochemical Profile

A biochemical profile was obtained for *E. coli-Xen14* using the api 20E system available from bioMérieux.

Sugar Utilization		Other Tests	
Glucose	+	α-galactosidase	+
Mannitol	+	Arginine Dihydrolase	-
Inositol	-	Lysine Decarboxylase	-
Sorbitol	-	Ornithine Decarboxylase	-
Rhamnose	+	Citrate Utilization	-
Sucrose	+	H <sub>2</sub> 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	Carbenicillin 100
Tetracycline 30	Kanamycin 30
Trimethoprim/ Sulfamethoxazole	Streptomycin 10
Gentamicin 10	Cephalothin 30
	Ampicillin 10

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

119223-R REV01