

Pseudomonas aeruginosa, P. aeruginosa PAO1 (Xen41)

Product Number: 119229

Material Provided: 1 Agar Plate

Storage Conditions: -80°C

Genetic Characteristics

Pseudomonas aeruginosa-Xen41 was derived from the parental strain *Ps. aeruginosa* PAO1, obtained from Pseudomonas Genetic Stock Center (PGSC). *P. aeruginosa* Xen41 possesses a single stable copy of the *P. luminescens luxCDABE* operon on the bacterial chromosome.

Growth Characteristics

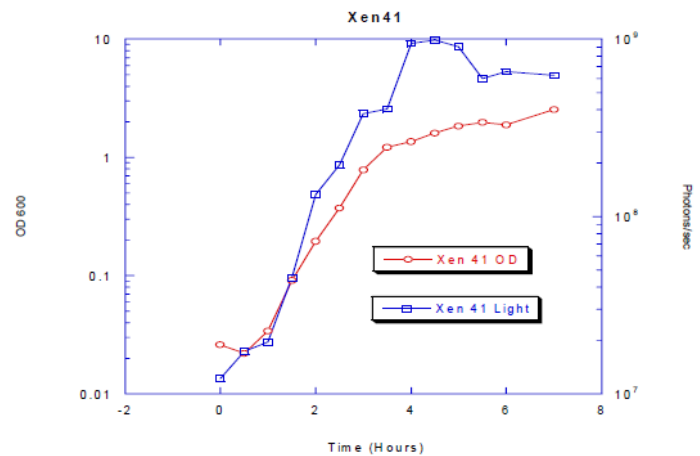
P. aeruginosa-Xen41 grows well in Luria Bertani (LB) medium at 37°C under ambient aeration. *P. aeruginosa*-Xen41 may also be grown selectively on LB agar containing 60 µg/mL Tetracycline.

Colonial Morphology

On LB plates, *P. aeruginosa*-Xen41 appears as large (3-5mm), yellow-green, irregularly round, mucoid colonies with butyrous centers after 24 hours incubation at 37°C.

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 3×10^8 cfu/mL of *Ps. Aeruginosa*-Xen41 and the relative light intensity is 0.013 photons/sec/cell.



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

119229-R REV01