

# APPLICATION NOTE

## VIBRIO CAMPBELLII QUORUM SENSING

### ABSTRACT

This study will demonstrate the use of the ATCC *Vibrio campbellii* Panel (ATCC® [MP-6™](#)) as a non-pathogenic model for AI-2-based quorum sensing pathways.

### INTRODUCTION

In many prokaryotes, cooperative behaviors are regulated through a density-dependent, signal-mediated communication system termed quorum sensing (QS).<sup>1</sup> When a bacterial population reaches a critical threshold, autoinducer signaling molecules (AI) specifically bind to a cognate regulatory protein or activate a two-component signal transduction system, leading to the regulation of group behaviors. In the marine organism *Vibrio campbellii*, AI signals (AI-1 and AI-2) and cognate regulators are used to regulate bioluminescence<sup>1</sup> (Figure 1). Since its discovery, AI-2 has proven ubiquitous within inter- and intraspecies communication, including that of pathogenic microorganisms.<sup>2</sup> Here, we show a panel of nine *V. campbellii* strains displaying wild-type or varying mutational phenotypes for use as a non-pathogenic model in the analysis of AI-2-based QS systems.

### MATERIALS AND METHODS

Nine *V. campbellii* strains were phenotypically analyzed for QS proficiency by monitoring the bioluminescence production of genotypically diverse strains that were plated together in pairs on Autoinducer Bioassay Medium.<sup>1,3-6</sup>

### RESULTS AND DISCUSSION

Upon analysis of paired strains, it was determined that bioluminescence could be restored in strains lacking regulator and/or AI production if the adjacent strain was proficient in that characteristic (Figure 2A-C, Table 1). Bioluminescence could not be restored in strains lacking part of the luxCDABE operon, which encodes for bioluminescence (Figure 1, Figure 2D, Table 1).

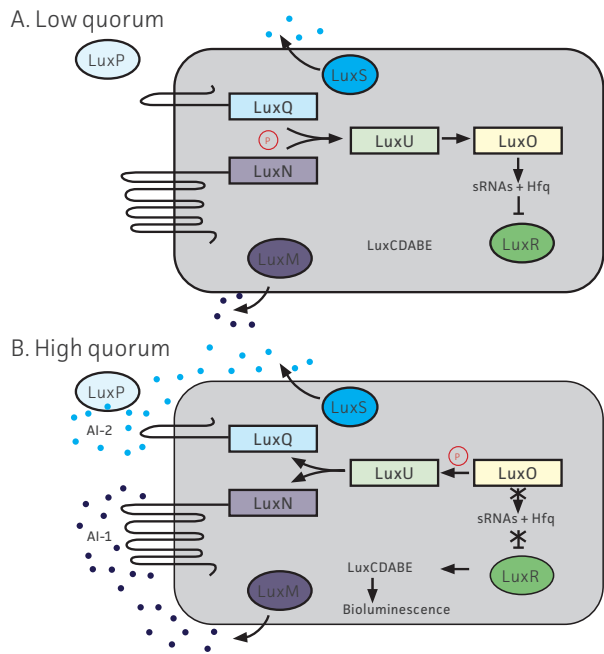


Figure 1: Quorum sensing

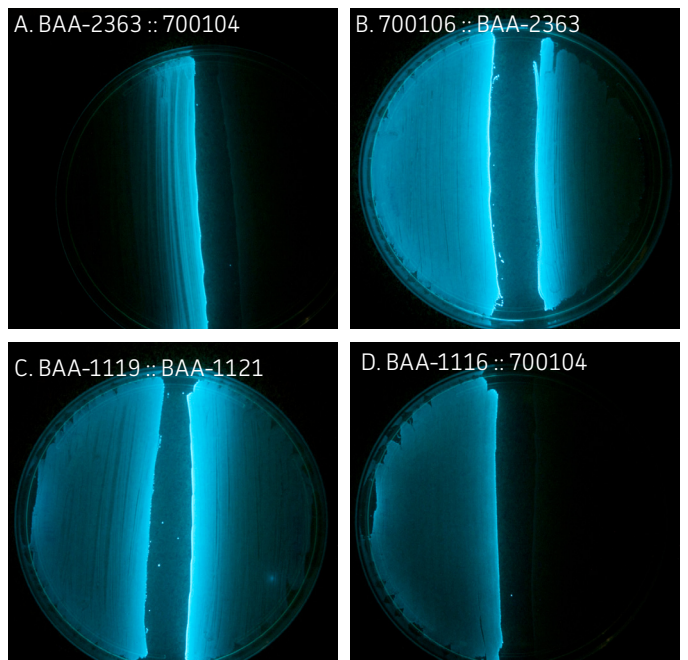


Figure 2: Bioluminescence

Table 1: ATCC *Vibrio campbellii* Panel (MP-6)

Sensors	ATCC No.	Autoinducers								
		1+, 2+	1+, 2+	1+, 2+	1+, 2+	1+, 2+	1-, 2+	1+, 2-	1+, 2-	1-, 2-
luxA-	700104™	-	-	-	-	-	-	-	-	-
1+, 2-	700106™	+	+	+	+	+	+	+	+	+
1+, 2+	BAA-1116™	+	+	+	+	+	+	+	+	+
1-, 2+	BAA-1117™	+	+	+	+	+	+	+	+	+
1+, 2-	BAA-1118™	+	+	+	+	+	+	+	+	+
1-, 2+	BAA-1119™	+	+	+	+	+	+	+	+	+
1+, 2+	BAA-1120™	+	+	+	+	+	+	+	+	+
1-, 2+	BAA-1121™	+	+	+	+	+	+	-	-	-
1+, 2+	BAA-2363™	+	+	+	+	+	-	-	-	-

Sensor 1 = LuxN; Sensor 2 = LuxQ; Autoinducer 1 = AI-1; Autoinducer 1 AI-2; (+) = Light observed; (-) = No Light observed


## CONCLUSION

The characterization of these *V. campbellii* strains illustrates that ATCC [MP-6](#) is well suited as a non-pathogenic model for the analysis of AI-2-based, two-component regulatory QS pathways.

## REFERENCES

- 1 Bassler B, Wright M, Silverman M. Multiple signaling systems controlling expression of luminescence in *Vibrio harveyi*: sequence and function of genes encoding a second sensory pathway. *Mol Microbiol* 13: 273-286, 1994.
- 2 Galloway WR, et al. Quorum sensing in Gram-negative bacteria: small-molecule modulation of AHL and AI-2 quorum sensing pathways. *Chem Rev* 111: 28-67, 2011.
- 3 Surette MG, Miller MB, Bassler BL. Quorum sensing in *Escherichia coli*, *Salmonella typhimurium*, and *Vibrio harveyi*: a new family of genes responsible for autoinducer production. *Proc Natl Acad Sci U S A* 96: 1639-1644, 1999.
- 4 Waters CM, Bassler BL. The *Vibrio harveyi* quorum-sensing system uses shared regulatory components to discriminate between multiple autoinducers. *Genes Dev* 20: 2754-2767, 2006.
- 5 Bassler BL, Greenberg EP, Stevens AM. Cross-species induction of luminescence in the quorum-sensing bacterium *Vibrio harveyi*. *J Bacteriol* 179: 4043-4045, 1997.
- 6 Bassler BL, Wright M, Silverman MR. Sequence and function of LuxO, a negative regulator of luminescence in *Vibrio harveyi*. *Mol Microbiol* 12: 403-412, 1994.
- 7 Lin B, et al. Comparative genomic analyses identify the *Vibrio harveyi* genome sequenced strains BAA-1116 and HY01 as *Vibrio campbellii*. *Environ Microbiol Rep* 2(1): 81-89, 2010.

 10801 University Boulevard  
Manassas, Virginia 20110-2209

 703.365.2700

 703.365.2701

 sales@atcc.org

 www.atcc.org

AP-122021-v05

©2022 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise. *Vibrio harveyi* ATCC BAA-1116 and its derived mutants have been reclassified as *Vibrio campbellii* based on microarray comparative genome hybridization and multilocus sequence analysis.

These products are for laboratory use only. Not for human or diagnostic use. ATCC products may not be resold, modified for resale, used to provide commercial services or to manufacture commercial products without prior ATCC written approval.