

SEEING IS BELIEVING - REPORTER LABELED MICROBIAL CONTROL STRAINS

Jodie Y. Lee, M.S.
Lead Biologist, ATCC
May 21, 2015

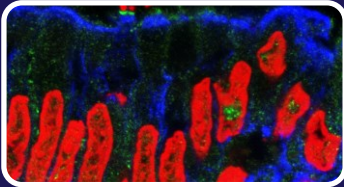


THE ESSENTIALS OF LIFE SCIENCE RESEARCH
GLOBALLY DELIVERED™

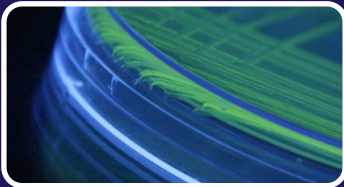
Today's discussion



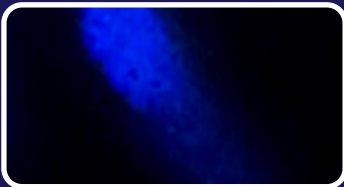
About ATCC



Reporter labels and their use



GFP-labeled strains



NanoLuc[®]-labeled Shiga toxin-producing *Escherichia coli* strains

About ATCC

- Founded in 1925, ATCC is a non-profit organization with headquarters in Manassas, VA
- World's premiere biological materials resource and standards development organization
- ATCC collaborates with and supports the scientific community with industry-standard products and innovative solutions
- Strong team of 400+ employees; over one third with advanced degrees
- Broad range of biomaterials
 - Cell lines, iPSCs, primary cells, and hTERT immortalized cells
 - Bacteria, yeasts, protists, and viruses
 - Tumor cell panels
 - Media, sera, and reagents

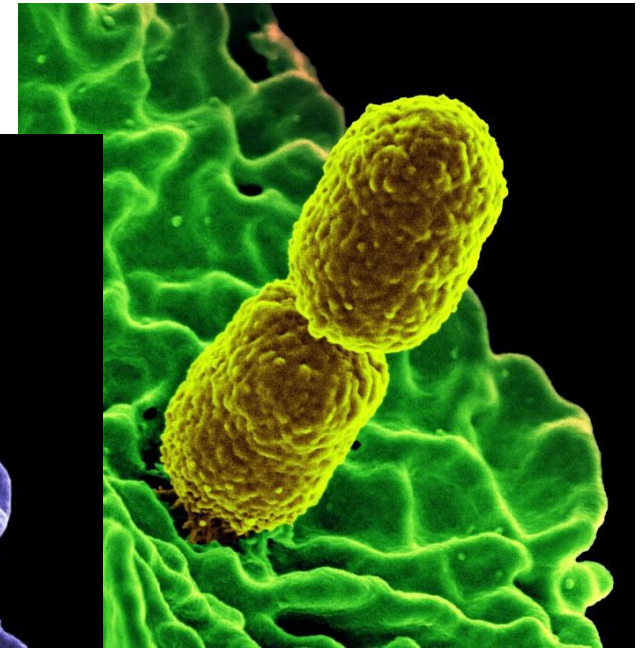
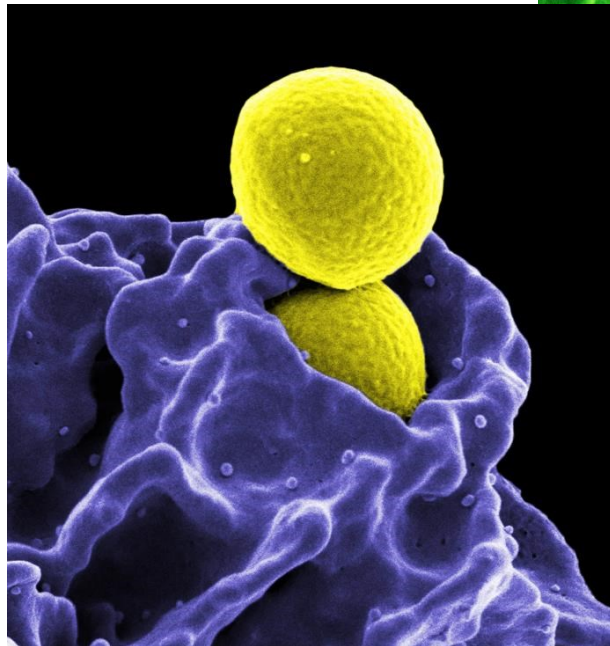


Established partner to global researchers and scientists

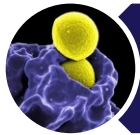


Why label bacteria?

- Visualize bacteria under experimental conditions
- Differentiate between sample contamination and false positives from control cross-contamination
 - Saves money, time, and worry



Why label bacteria?



Microbial quantification and detection



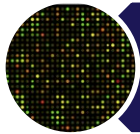
Host-pathogen interactions



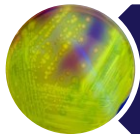
Drug discovery



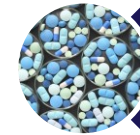
Food testing



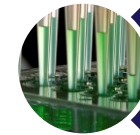
Gene expression



Biochemical interactions

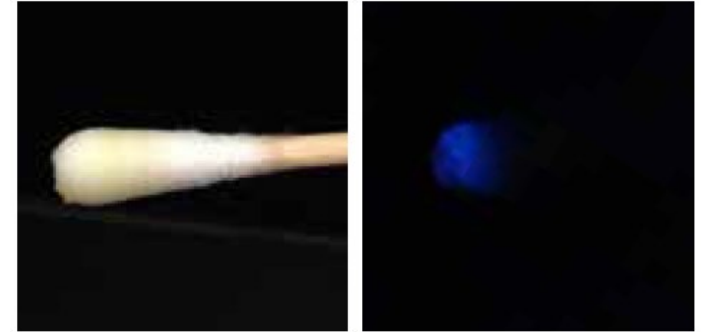
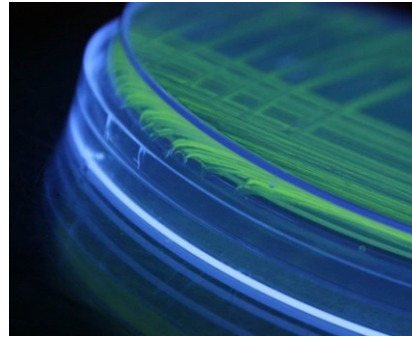


Novel therapeutics



Quality control assays

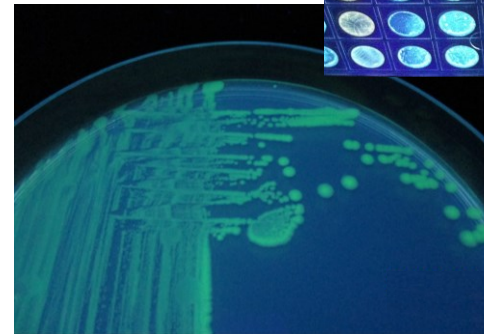
Reporter characteristics



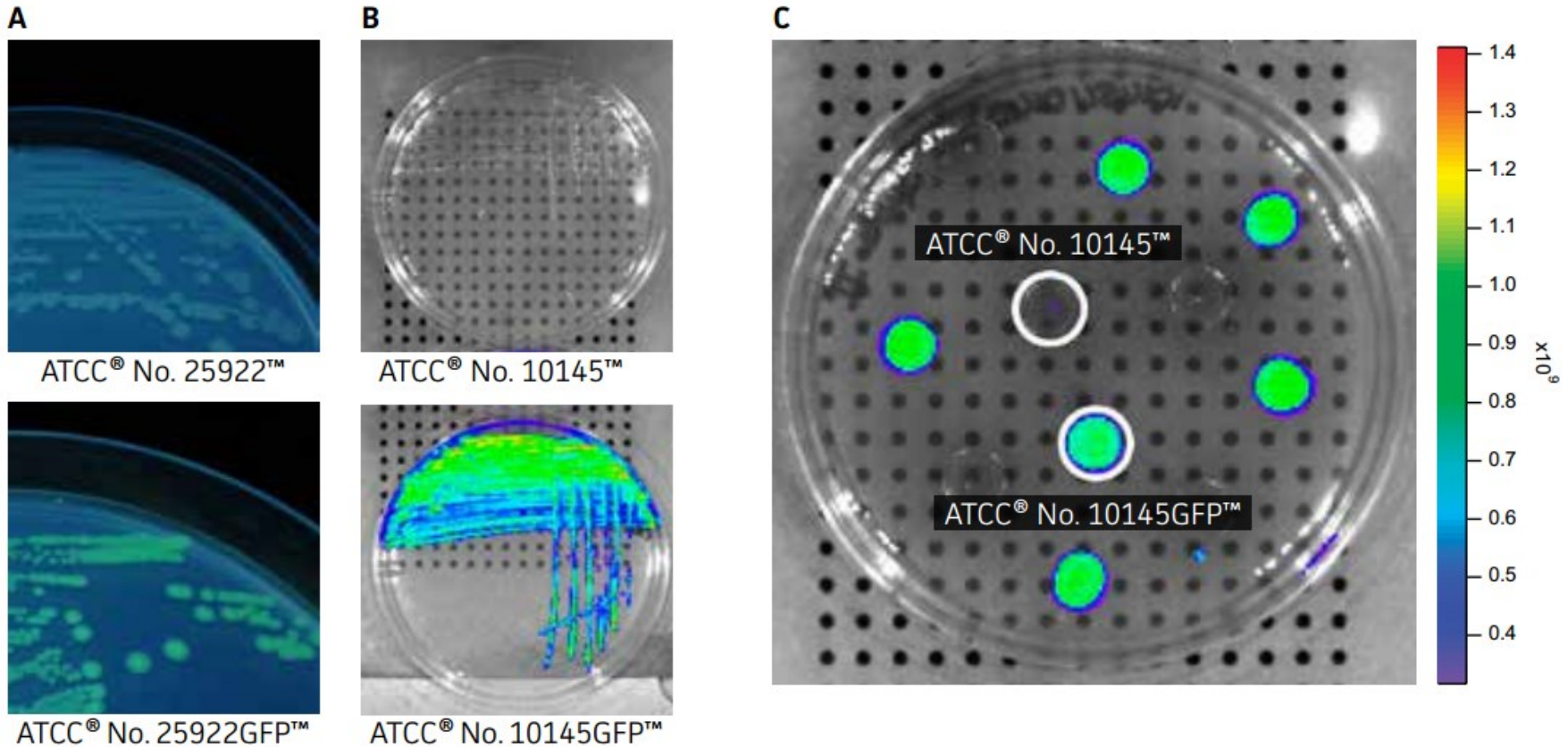
	Green Fluorescent Protein (GFP)	NanoLuc [®] Luciferase
Live-cell assays	Yes	No
Well-characterized	Yes	No
Highly stable	Yes	Yes
Substrate required	No	Yes
Glow response	Yes	Yes
Use <i>in vivo</i>	Yes	No

GFP-labeled strains

- Labeled Gram-negative organisms with GFPmut3a
 - *Escherichia coli*
 - *Salmonella enterica* subsp. *enterica* serovar Typhimurium
 - *Shigella flexneri*
 - *Pseudomonas aeruginosa*
- *gfp* expressed behind the P_{lac} constitutive promoter
- Bright, consistent label can be used for:
 - Microbial quantification and detection
 - Host-pathogen interactions
 - Drug discovery
 - Food testing

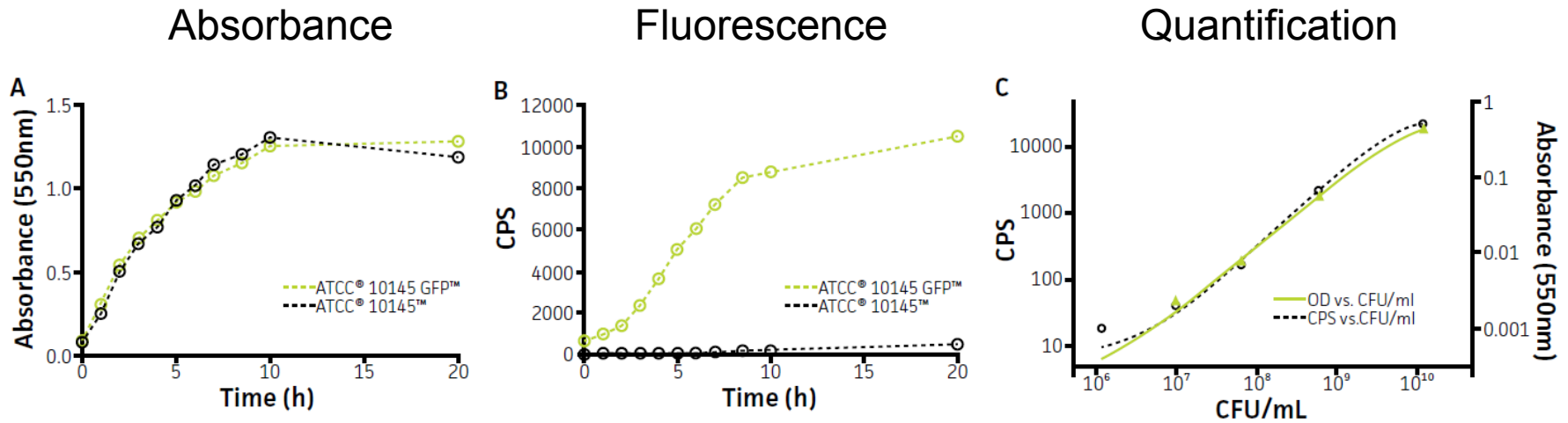


Visual detection of GFP-labeled organisms



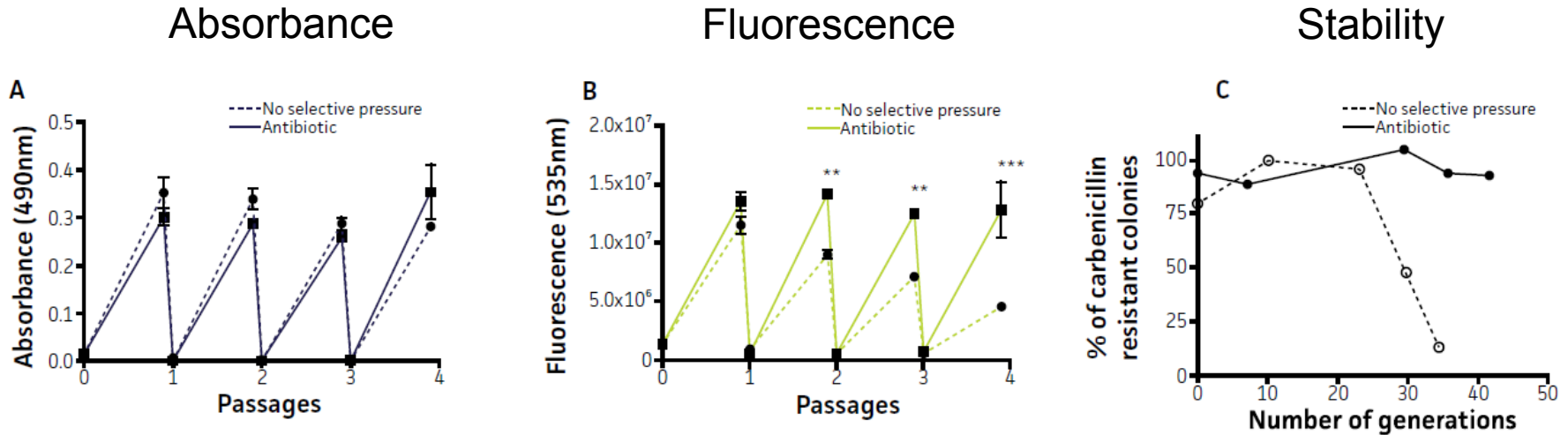
The expression of a bright GFP variant on a high-copy number plasmid facilitates visual identification when exposed to UV light (A) or imaged using a detection system (B & C)

Bacterial fitness and quantification



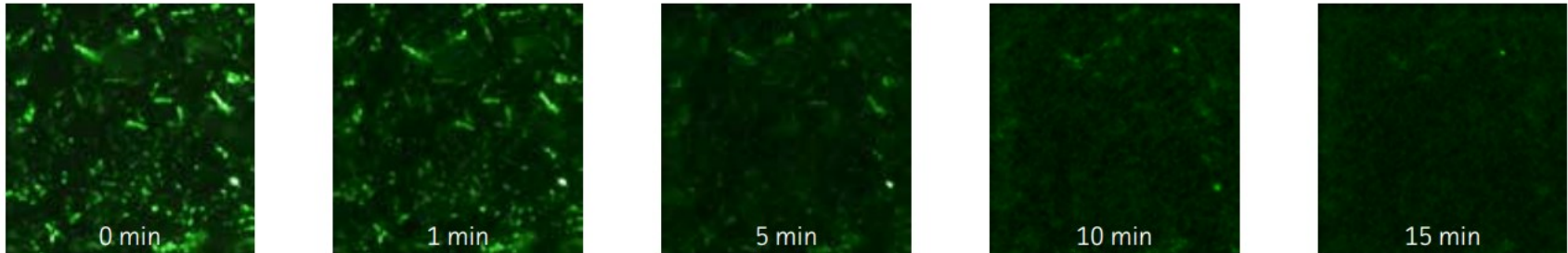
- GFPmut3a is not detrimental to bacterial fitness
- Fluorescence can be used for bacterial quantification

Plasmid stability

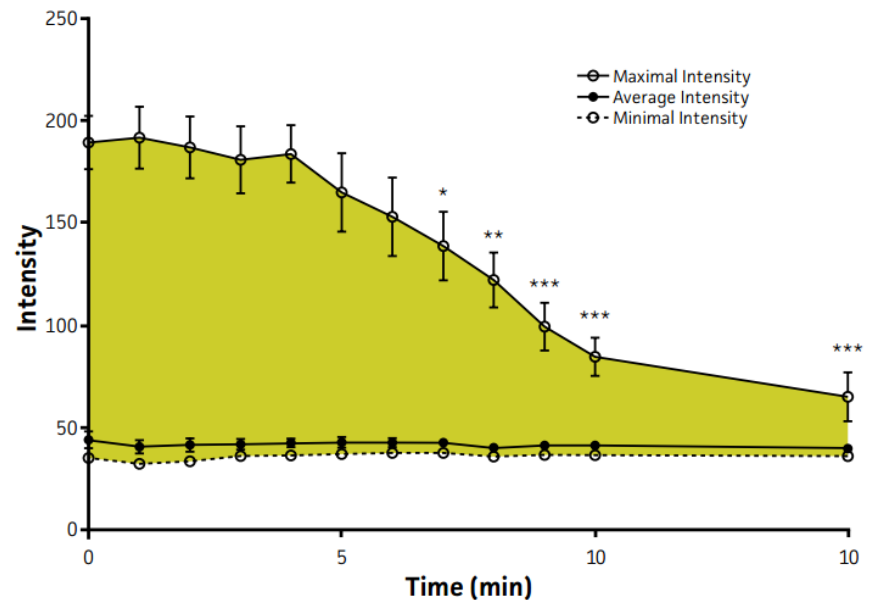


Plasmid is stable for at least 20 generations at 37°C in the absence of antibiotic pressure

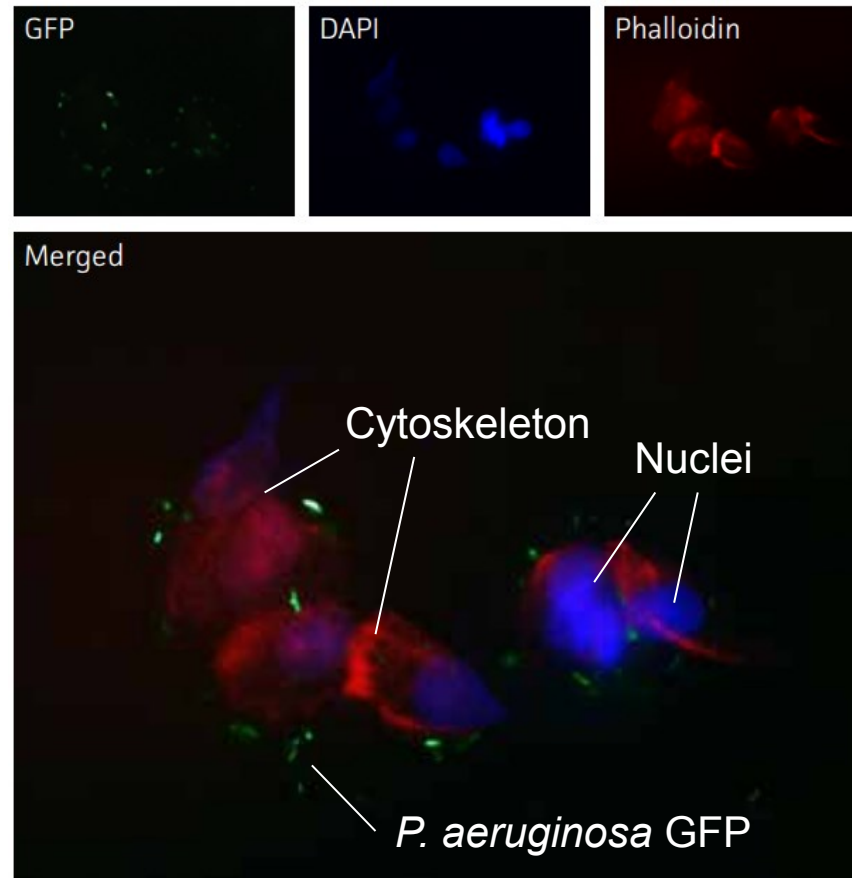
Photo bleaching



- *P. aeruginosa*-GFP (ATCC® 10145GFP™) was continuously exposed to UV light to determine resistance to photo bleaching
- Fluorescence remained stable during the first 5 minutes
- Significant loss of signal after 7 minutes of exposure



Fluorescent microscopy detection of pathogen-host interactions

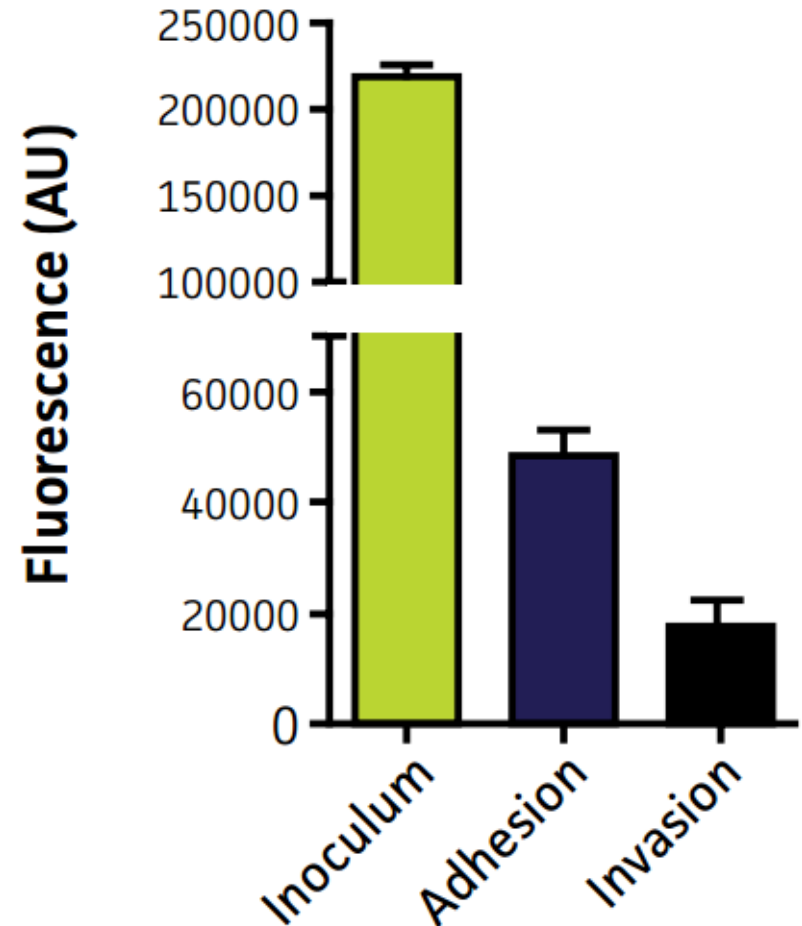


P. aeruginosa GFP (ATCC®10145GFP™) interaction with A549 (ATCC® CCL-185™) airway epithelial cells (100X magnification)

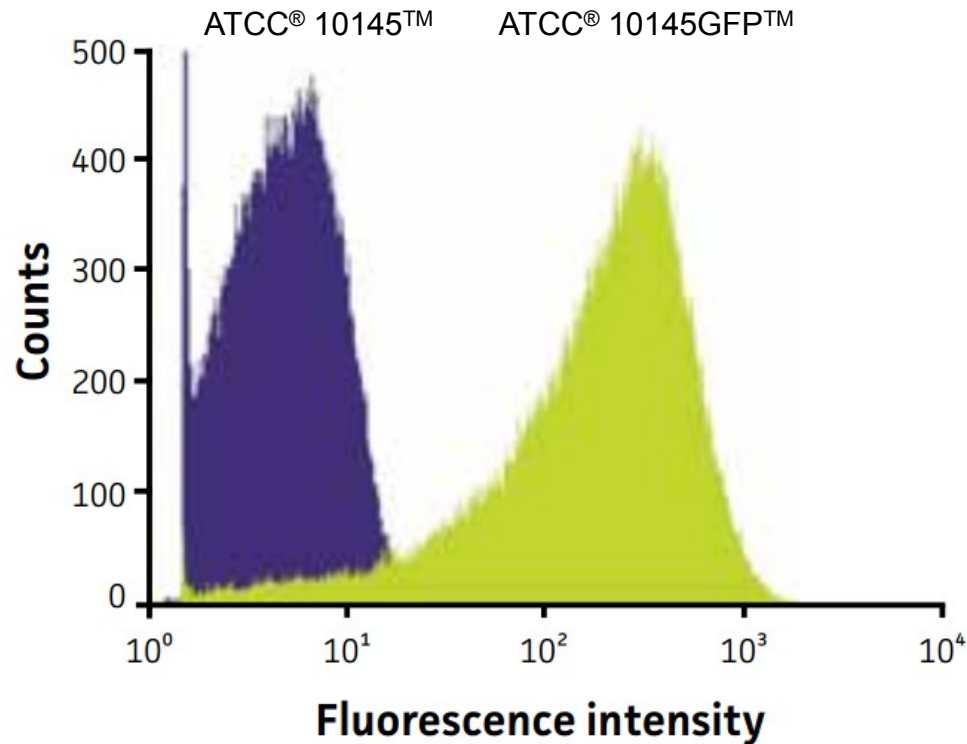
High-throughput detection of pathogen-host interactions

Invasion study

- *P. aeruginosa* GFP (ATCC®10145GFP™) was incubated in the presence of A549 (ATCC® CCL-185™) airway epithelial cell monolayers
- Cells were washed and medium supplemented with 100 µL gentamicin was added to kill extracellular bacteria (Adhesion)
- Cells were washed again (Invasion) and measured on a microplate reader



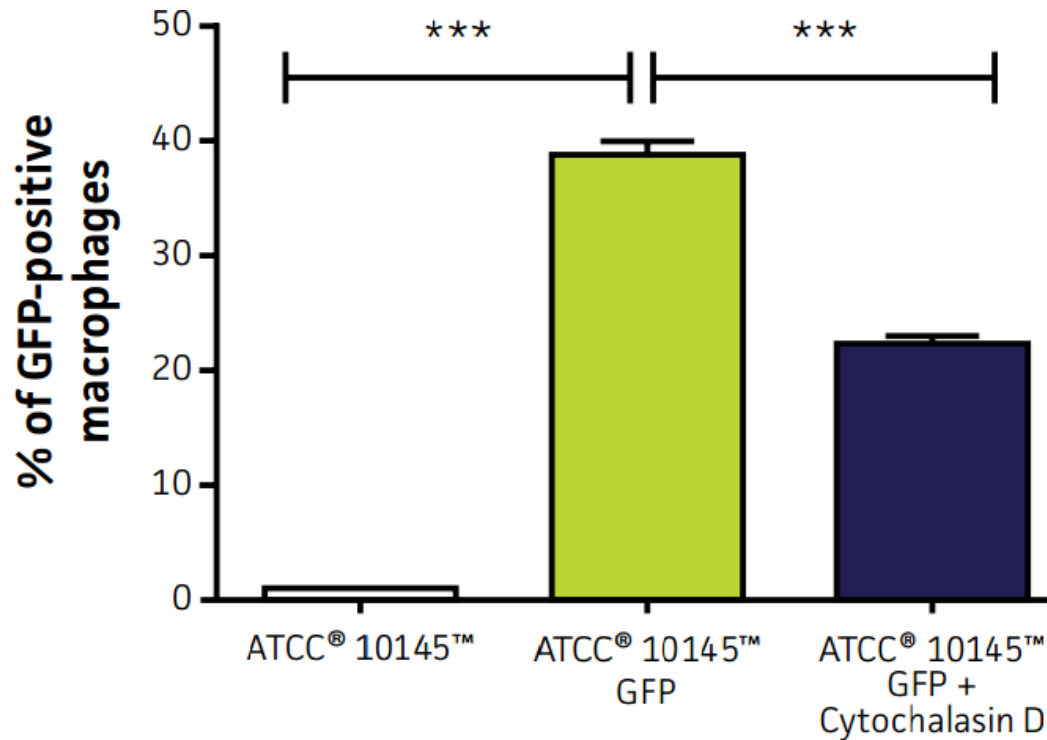
High-throughput detection of pathogen-host interactions



Flow Cytometry

- *P. aeruginosa* (ATCC® 10145™) (purple) and *P. aeruginosa* GFP (ATCC® 10145GFP™) (green) suspensions were analyzed by flow cytometry

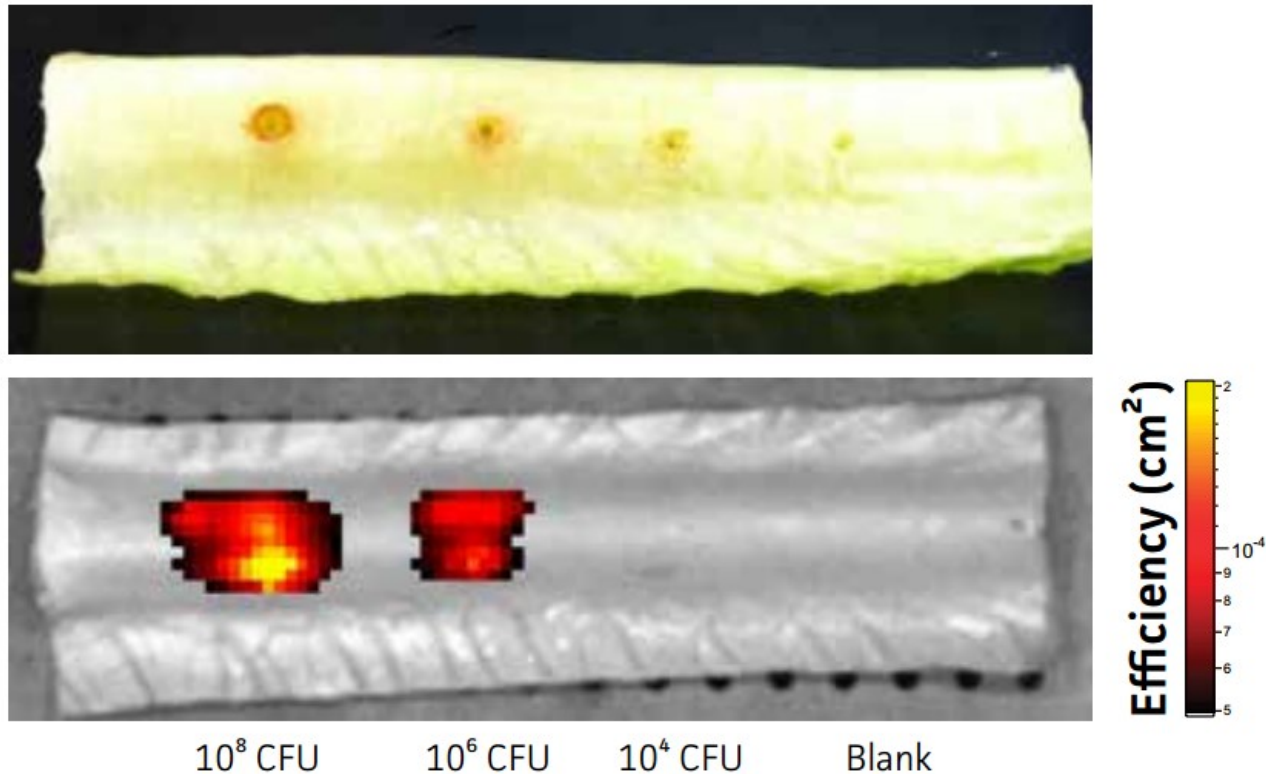
High-throughput detection of pathogen-host interactions



Macrophage uptake of *Pseudomonas aeruginosa*

- *P. aeruginosa* (ATCC® 10145™) and *P. aeruginosa* GFP (ATCC® 10145GFP™) were incubated in the presence or absence of Cytochalasin D and analyzed by flow cytometry

In vivo detection of GFP-labeled *P. aeruginosa*



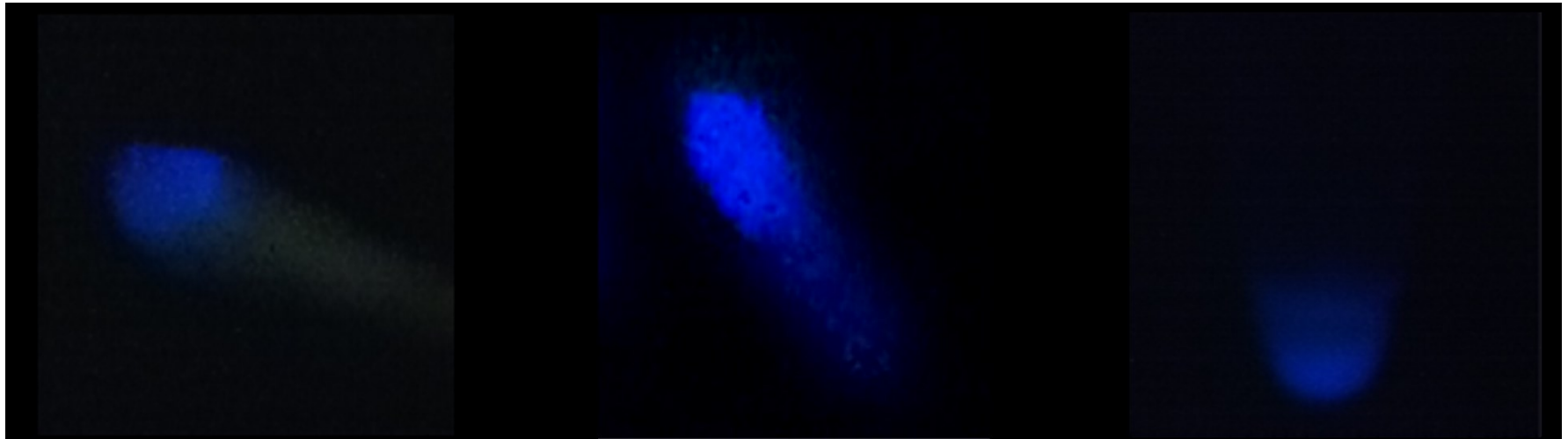
- Various doses of *P. aeruginosa* GFP (ATCC[®] 10145GFP[™]) were injected into the mid-rib of *Lactuca sativa* L. var. *longifolia*
- The bacteria was easily detected at higher concentrations in the plant host, indicating that this vector can be successfully employed to monitor bacterial growth within a plant host

ATCC GFP-labeled strains

ATCC® No.	Species	Reporter	Parental Strain (ATCC® No.)
25922GFP™	<i>Escherichia coli</i>	GFP	25922™
14028GFP™	<i>Salmonella enterica</i> subsp. <i>enterica</i> serovar Typhimurium	GFP	14028™
12022GFP™	<i>Shigella flexneri</i>	GFP	12022™
10145GFP™	<i>Pseudomonas aeruginosa</i>	GFP	10145™
15692GFP™	<i>Pseudomonas aeruginosa</i>	GFP	15692™

www.atcc.org/reporters

Promega NanoLuc[®] reporter



Cotton Swab

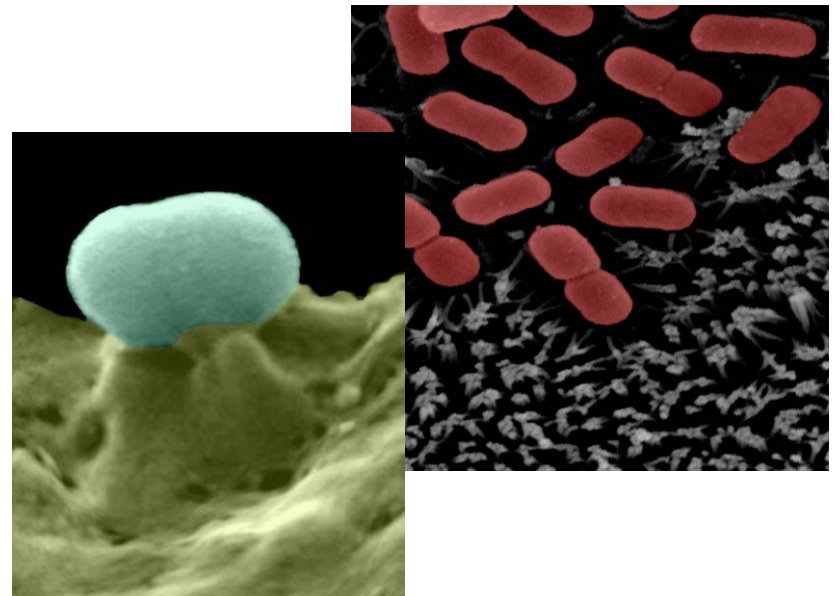
Filter Paper

Cell Pellet

- Intensely bright reporter
- Glow response
- Portable, does not require instrumentation

Shiga toxin-producing *E. coli*

- >265,000 cases of STEC infection in the United States each year
 - *E. coli* O157:H7 accounts for about 36% of STEC infections
 - ~5-10% of diagnosed infections develop into hemolytic uremic syndrome, a life threatening complication which can cause permanent health damage
- Food Safety and Modernization Act calls for expanded testing to include Non-O157 strains
 - O26
 - O45
 - O103
 - O111
 - O121
 - O145



Microbial strain authentication

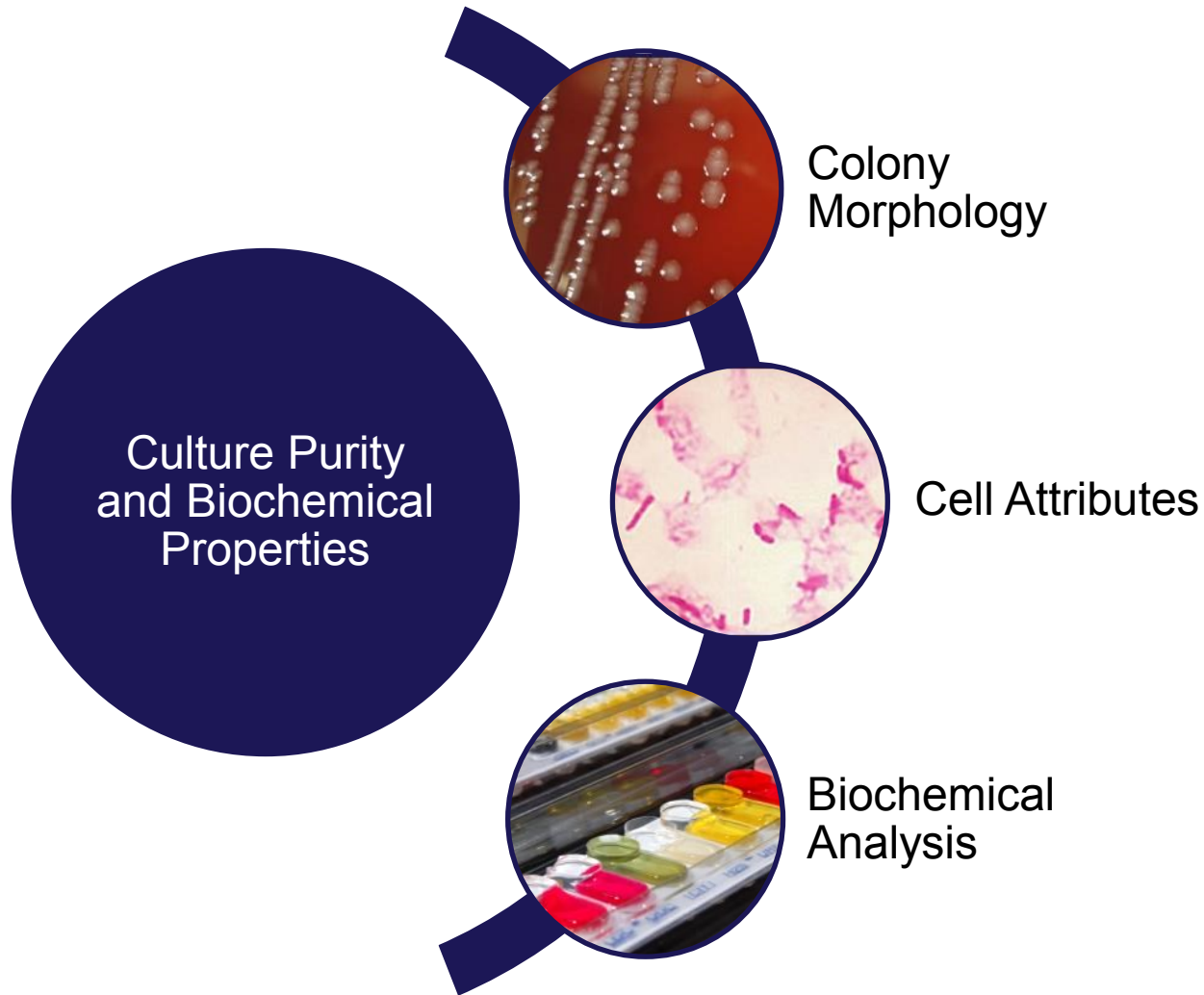


ATCC utilizes both classical and modern techniques

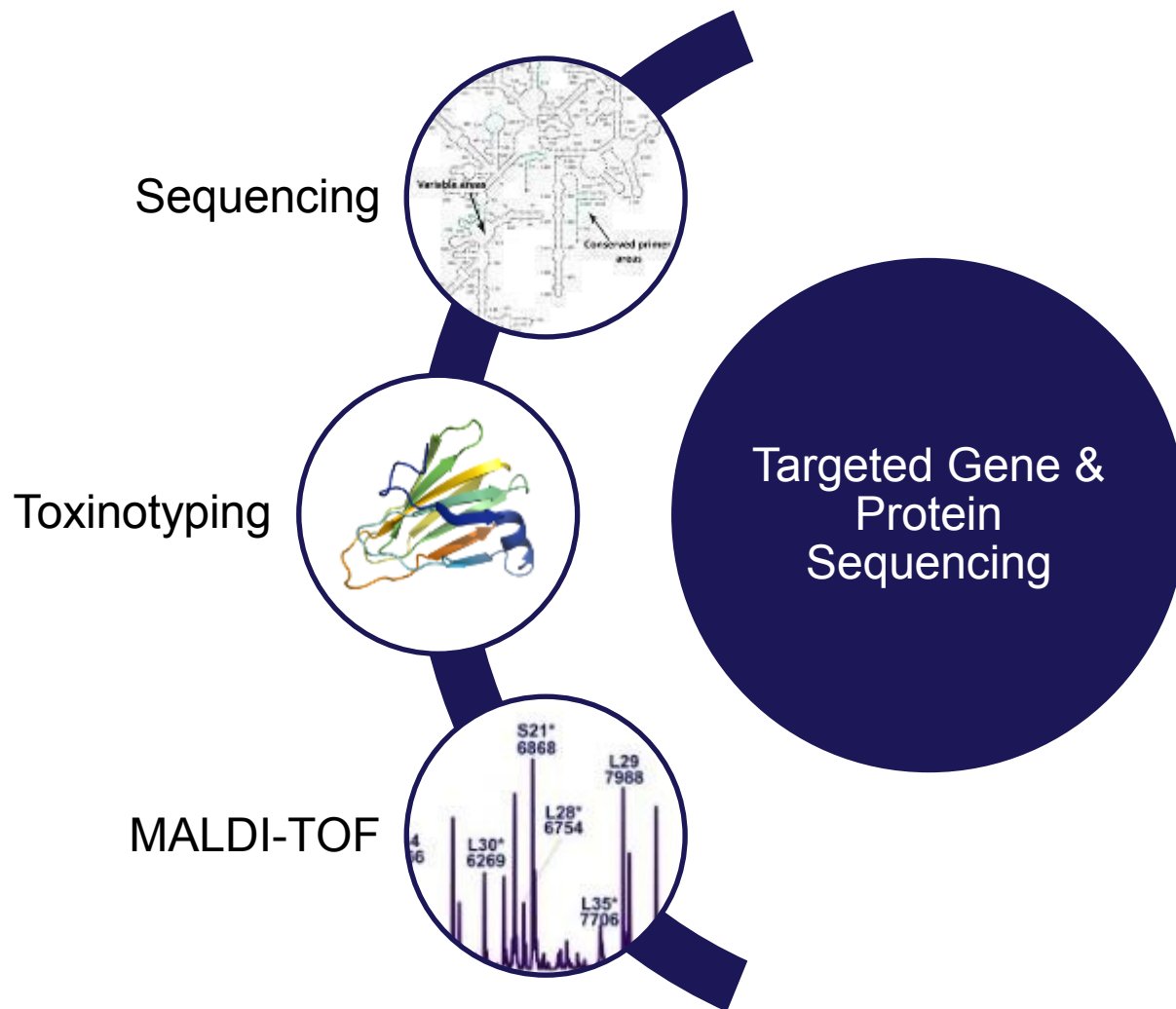
- Phenotypic analysis
- Genotypic & proteotypic analyses
- Functional analysis

No single method of identification is sufficient

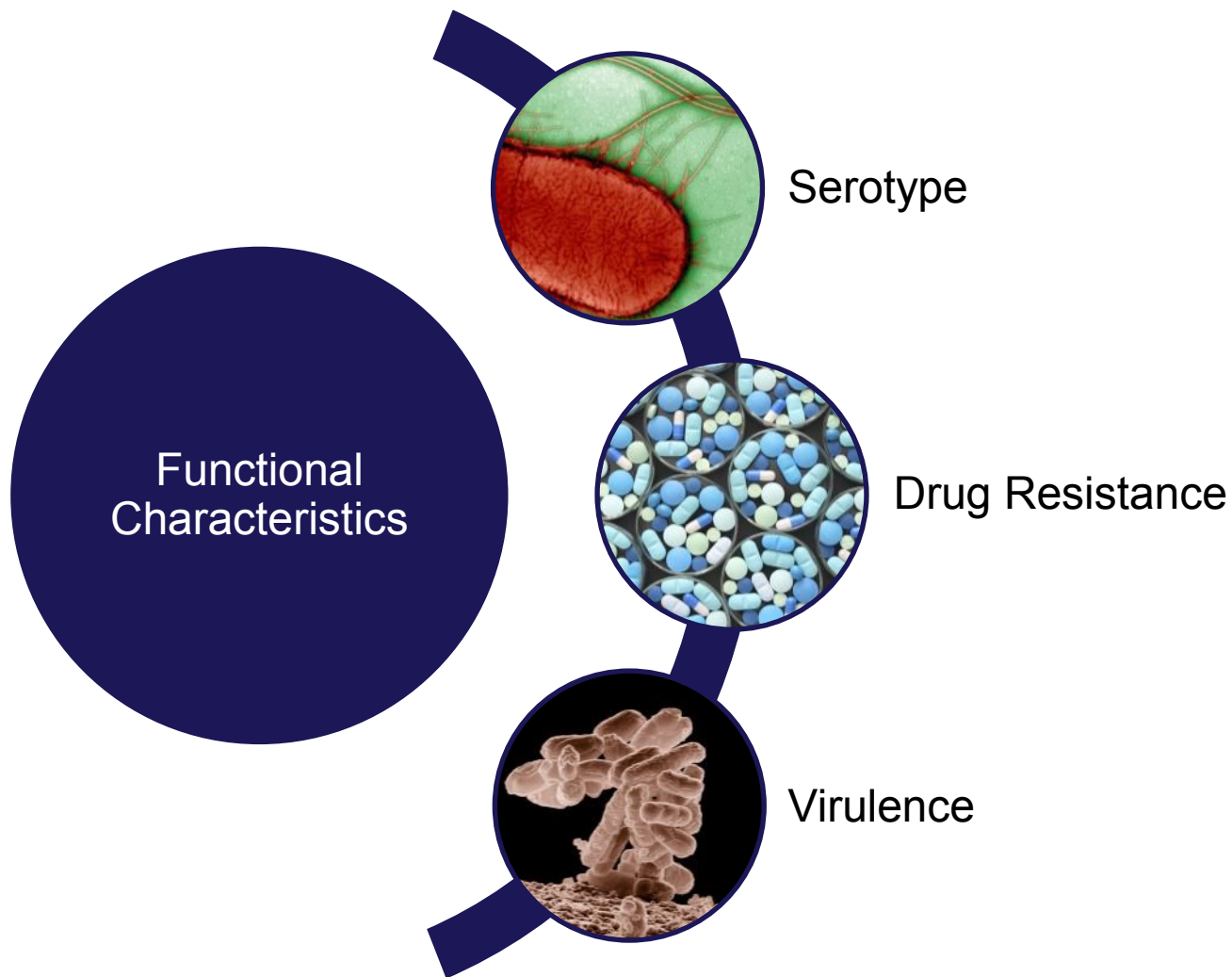
Phenotypic testing



Genotypic & proteotypic testing



Functional testing

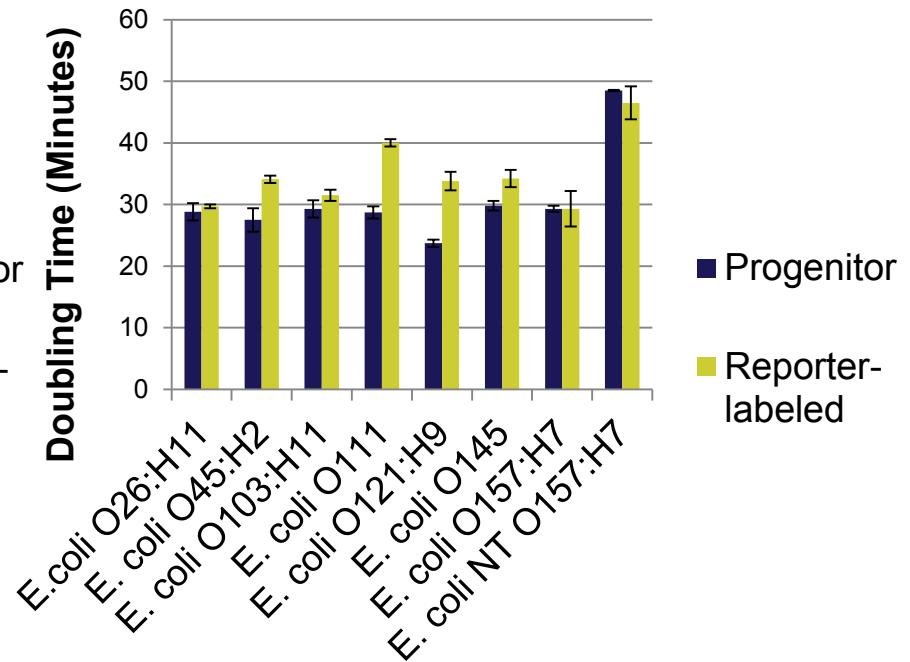
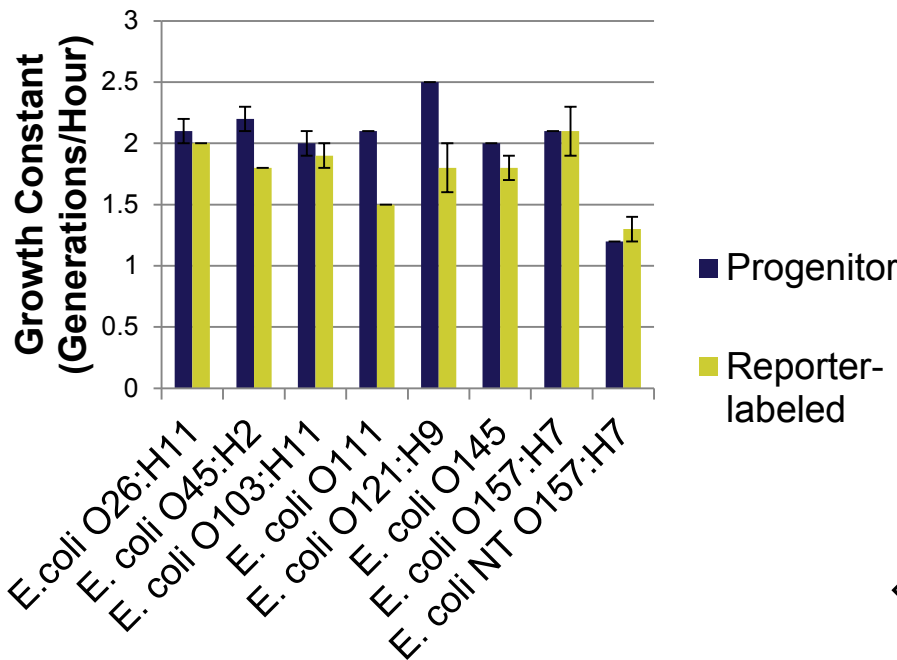


STEC Testing

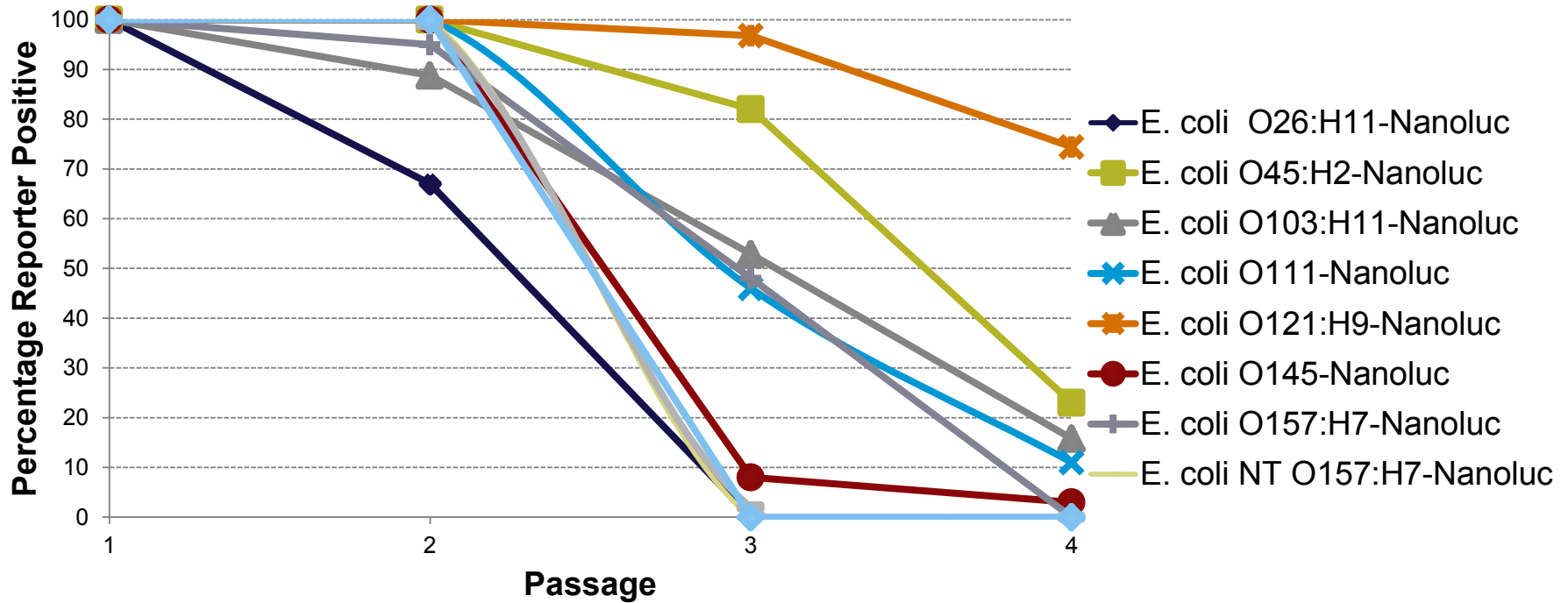
- Bacterial identification
 - VITEK[®] 2, VITEK[®] MS, API[®] Strips, 16S rRNA sequencing
- Molecular characterization to assess the presence of the *stx1*, *stx2*, and *eaeA* genes
 - PCR
- Serogroup identification
 - Immunoprecipitation assay



Minimal reporter effects on bacterial fitness



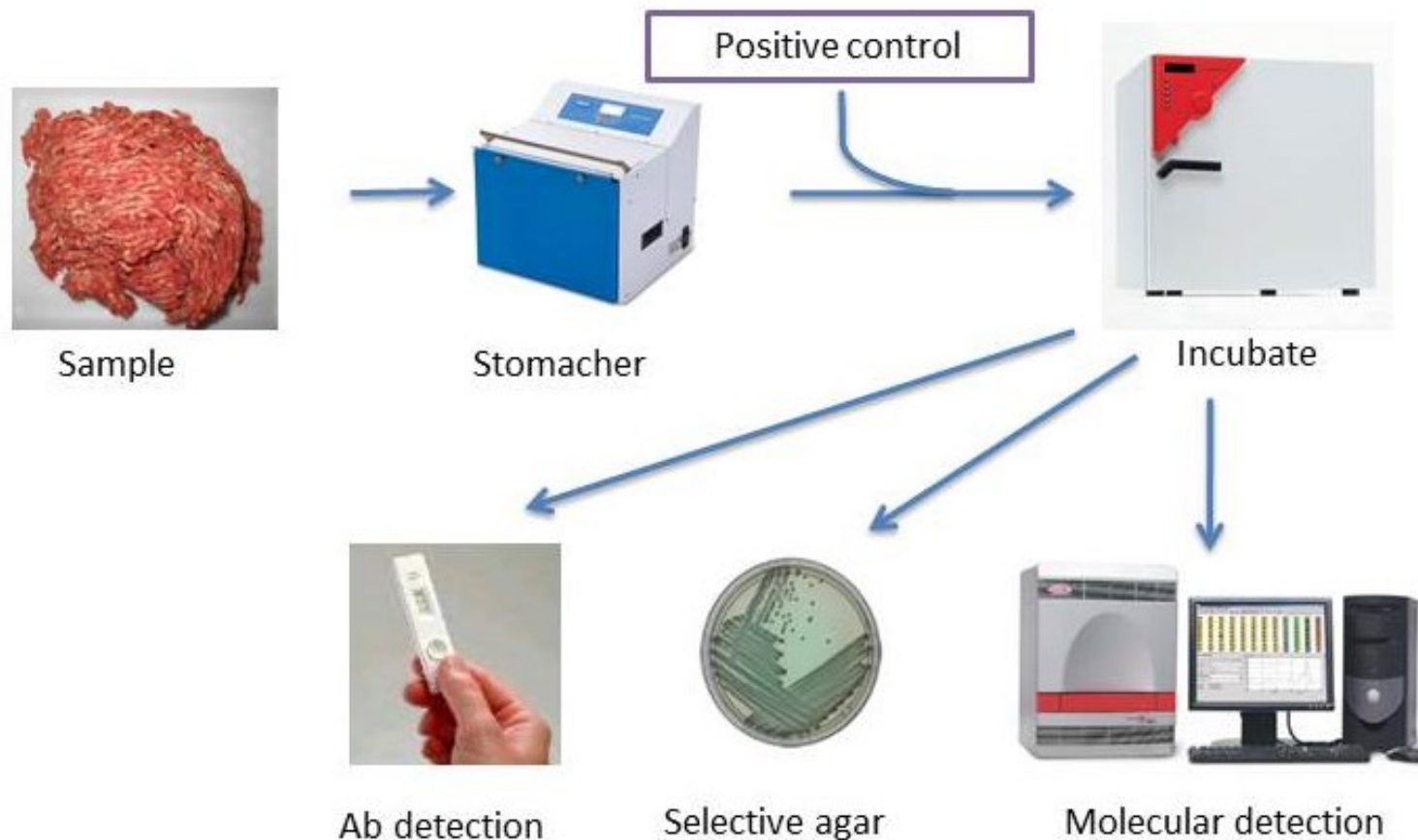
Plasmid stable for ≥ 2 days



Observe reporter in $>65\%$ of c.f.u. after 2 days at 42°C

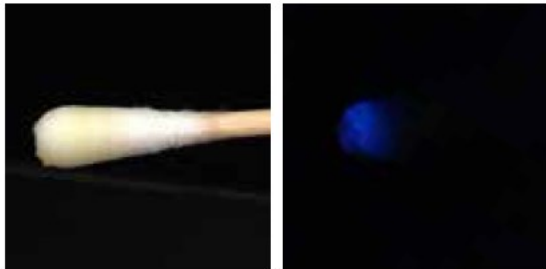
Shiga toxin-producing *E. coli* (STEC) control strains

Labeled controls rule out cross-contamination of samples with the control strain

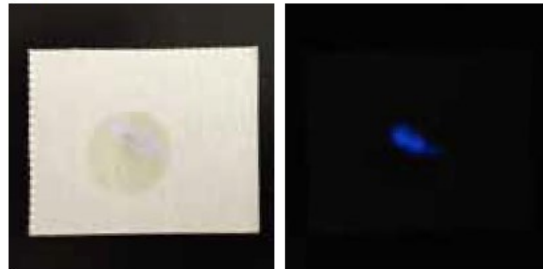


NanoLuc[®]-labeled control detection

Cotton Swab



Paper



Broth



External testing confirmed a clear signal after FSIS protocol for ground beef

NanoLuc[®]-labeled *Escherichia coli*

ATCC [®] No.	Species	Serotype	Genotype
BAA-2580-PACK [™]	<i>Escherichia coli</i>	O26:H11	<i>stx1+</i> , <i>stx2+</i> , <i>eaeA+</i>
BAA-2581-PACK [™]	<i>Escherichia coli</i>	O45:H2	<i>stx1+</i> , <i>stx2-</i> , <i>eaeA+</i>
BAA-2582-PACK [™]	<i>Escherichia coli</i>	O103:H11	<i>stx1+</i> , <i>stx2-</i> , <i>eaeA+</i>
BAA-2583-PACK [™]	<i>Escherichia coli</i>	O111	<i>stx1+</i> , <i>stx2-</i> , <i>eaeA+</i>
BAA-2584-PACK [™]	<i>Escherichia coli</i>	O121:H19	<i>stx1-</i> , <i>stx2+</i> , <i>eaeA+</i>
BAA-2585-PACK [™]	<i>Escherichia coli</i>	O145	<i>stx1-</i> , <i>stx2+</i> , <i>eaeA+</i>
BAA-2586-PACK [™]	<i>Escherichia coli</i>	O157:H7	<i>stx1+</i> , <i>stx2+</i> , <i>eaeA+</i>
BAA-2587-PACK [™]	<i>Escherichia coli</i>	O157:H7	<i>stx1-</i> , <i>stx2-</i> , <i>eaeA-</i>

www.atcc.org/reporters

Coming soon – Big-Six *Escherichia coli* NanoLuc[®] Strains Panel

Additional testing resources



Food Testing
Reference
Strains

www.atcc.org/Food



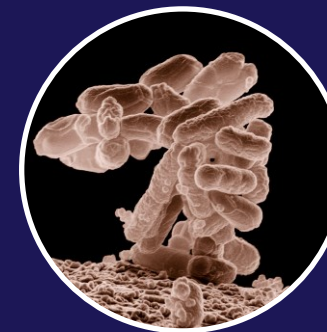
Proficiency
Testing
Programs

www.atcc.org/PTPrograms



ATCC® Minis
Quality Control
Strains

www.atcc.org/Minis



“Big Six” STEC
DNA and Strains
Panels

www.atcc.org/MP

ATCC Food Testing Solutions



Conclusions

- Reporter labels are a flexible research tool
- Using labels for microbial detection enhances safety and reliability in food testing
- ATCC strains provide
 - Authenticated reference standards
 - Low Passage
 - Significant savings in time and effort

Acknowledgements

- Mariette Barbier, Ph.D.

© 2015 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. NanoLuc[®] is a registered trademark of Promega Corporation. VITEK[®] and API[®] are registered trademarks of bioMérieux.

Thank you!

Watch recorded ATCC® “*Excellence in Research*” webinars on demand at www.atcc.org/webinars.



The ATCC® “*Excellence in Research*” webinar series returns in Fall 2015. Look for webinars starting in September at www.atcc.org/webinars.

Thank you for joining today!
Please send additional questions to tech@atcc.org