



Verification Testing of iQ-Check™ and dd-Check STEC Kits with Onions, Carrots, and Cucumbers

Jennifer Pelowitz, Wendy Lauer, and Joshua Whitworth
Bio-Rad Laboratories, Inc., Hercules, CA 94547 USA

Abstract

While fresh produce is nutritionally beneficial, it can pose a risk of foodborne outbreaks due to bacterial contamination, especially because it is often consumed raw. Food safety standards and preventive measures are needed to protect the safety of these commodities. The purpose of this study was to perform verification testing on onions, carrots, and cucumbers with iQ-Check PCR Detection Kits and dd-Check STEC Kits (onions only) to determine their fit for purpose for these matrices.

Introduction

Salmonella spp. and *Escherichia coli* are two of the most common bacterial pathogens responsible for foodborne outbreaks. Both can contaminate a wide range of foods, leading to significant public health concerns. In general, they can be present in raw fruits and vegetables due to contamination during farming, harvesting, or processing or due to contaminated water sources.

In a survey of foodborne illness from 1998 to 2022, the Centers for Disease Control and Prevention (CDC) attributed 66.8% and 39.4% cases of *E. coli* and *Salmonella* illness, respectively, to a fruit or vegetable food matrix. Over the last few months of 2024, there were outbreaks in the United States associated with organic carrots, cucumbers, and onions, with the last being linked to a popular fast-food chain. These three outbreaks alone involved 265 reported cases and led to 82 hospitalizations.

The use of molecular methods for pathogen detection are useful for rapid determination of results, particularly for produce matrices that have a short shelf life. Real-time PCR (qPCR) has the added advantage of increased sensitivity for the target organism and selectivity against competing bacteria. Droplet Digital™ PCR (ddPCR™) technology is a powerful tool in which a sample is divided into thousands of nanoliter-sized droplets. This approach enables direct colocalization, which is useful for detection and confirmation of, for example, Shiga toxin–producing *E. coli* (STEC), ensuring the target virulence genes *stx1* and *stx2*, as well as *eae*, are detected as physically present on the genome of a single bacterial cell and each not merely present in different cells in the sample.

In this study, we performed verification testing to determine the viability of iQ-Check PCR Kits for detecting pathogens in onions, carrots, and cucumbers. Onion samples were also tested with the dd-Check STEC Kit.

Methods

Matrices

Fresh chopped onions, sliced carrots, and whole cucumbers were purchased from a local grocery store.

Test Strains

Cultures were prepared by inoculating 9 ml Tryptic Soy Broth (TSB; Hardy Diagnostics, catalog #88) and incubating for 18 hr at $37 \pm 1^\circ\text{C}$, followed by refrigeration at $4 \pm 1^\circ\text{C}$. Cultures were serially diluted and plated on Tryptic Soy Agar (TSA; Hardy Diagnostics, #G60) for 18 hr before inoculating matrices to determine titers for target inoculation. Cultures were plated on TSA again at the time of inoculation. Target strains were paired with competing organisms for inoculation (Table 1).

Enrichment Conditions and PCR/ddPCR Methods

The onion and carrot samples were enriched with Buffered Peptone Water Standard (BPW Standard; Bio-Rad Laboratories, Inc., #12013259) prewarmed to $37 \pm 1^\circ\text{C}$ at a 1:9 ratio (25 g sample + 225 ml medium). Selective STEC Supplement (Bio-Rad, #3564005) was added to half of the onion enrichments. Samples were incubated for 18 hr at $37 \pm 1^\circ\text{C}$. The iQ-Check Easy II DNA extraction protocol was used for samples tested with the iQ-Check *E. coli* O157:H7 PCR Detection Kit (Bio-Rad, #3578114) and iQ-Check STEC VirX PCR Detection Kit (Bio-Rad, #3578139).

Table 1. Target organisms matched with competing strains.

Food Matrix	Target Organism	Strain ID	Competing Organism	Strain ID
Onion	<i>E. coli</i> O157:H7	FDA ESC1177	<i>E. coli</i>	ATCC 25922
Carrot	<i>E. coli</i> O157:H7	FDA ESC1177	<i>E. coli</i>	ATCC 25922
Cucumber	<i>S. Typhimurium</i>	ATCC 14028	<i>E. coli</i>	ATCC 25922

The tube filtration and centrifugation protocols were used for samples tested with the dd-Check STEC Kit (Bio-Rad, #12009999). PCR was run on either the CFX96 Touch Deep Well or the CFX Opus Deepwell (Bio-Rad, #17007991) Real-Time PCR System. For Droplet Digital PCR, droplets were generated with the QX200™ Droplet Generator (Bio-Rad, #17005227) and read with the QX200 Droplet Reader (Bio-Rad, #17005228).

Whole cucumbers were sliced and enriched with BPW Standard prewarmed to 37 ± 1°C at a 1:9 ratio (25 g sample and 225 ml medium). Samples were incubated for 18 hr at 37 ± 1°C. The iQ-Check Easy I DNA extraction protocol was used, and samples were tested with the iQ-Check *Salmonella* II Kit (Bio-Rad, #3578123).

Test Protocol

Cultures were prepared based on the 18 hr–incubation plate count, and the final count was calculated by plating the refrigerated culture used for inoculation. For high level–inoculum samples, the targeted bacteria range was 100–1,000 colony forming units (CFU), and the competing bacteria range was >1,000 CFU. For low level–inoculum samples, the targeted bacteria range was 5–10 CFU, and the competing bacteria range was 100–500 CFU. Any deviation from these values was noted.

For each matrix tested, three samples were spiked with high-level inoculum of target bacteria and competing bacteria. In addition, three samples were spiked with low-level inoculum of target and competing bacteria; one sample not inoculated was used as a matrix control.

After 18 hr of enrichment, samples were removed from the incubator and tested with iQ-Check PCR and dd-Check STEC Kits.

Results

Detection of *E. coli* O157:H7 in Onions

When tested in triplicate with the iQ-Check *E. coli* O157:H7 PCR Kit, positive results were obtained for all inoculation levels after 18 hr of incubation, with or without STEC Supplement (Table 2). All inoculation levels also yielded positive results after 18 hr of incubation, with or without STEC Supplement, when tested with the iQ-Check STEC VirX PCR Kit (Table 3). When tested with dd-Check STEC Solution, all inoculation levels showed strong positive results after 18 hr of incubation. The samples were run with dithiothreitol (DTT; Bio-Rad, #12012171), and dilution from 1:100 to 1:1,000 was needed to ascertain linkage (Table 4, Figure 1).

Table 2. iQ-Check *E. coli* O157:H7 PCR Kit results for onions.

Matrix	Inoculum	Sample	Rep 1			Rep 2			Rep 3			PCR Curves		
			FAM	HEX	Result	FAM	HEX	Result	FAM	HEX	Result			
Onions, fresh chopped 18 hr incubation BPW	No inoculum	0	N/A	33.5	Negative	N/A	32.98	Negative	N/A	33.66	Negative			
	<i>E. coli</i> O157:H7: 560 CFU <i>E. coli</i> : 3,794 CFU	H1	18.03	34.25	Positive	17.85	34.55	Positive	17.88	34.45	Positive			
		H2	17.95	34.14	Positive	17.74	34.14	Positive	17.61	34.49	Positive			
		H3	17.51	34.39	Positive	17.48	35.96	Positive	17.46	34.87	Positive			
	<i>E. coli</i> O157:H7: 6 CFU <i>E. coli</i> : 379 CFU	L1	19.81	33.84	Positive	19.82	33.36	Positive	20.12	32.86	Positive			
		L2	22.80	31.68	Positive	22.94	31.85	Positive	23.02	32.50	Positive			
		L3	20.07	33.03	Positive	20.07	33.08	Positive	19.98	33.50	Positive			
	Onions, fresh chopped 18 hr incubation BPW + STEC Supplement	No inoculum	0	N/A	33.34	Negative	N/A	33.01	Negative	N/A	33.16		Negative	
		<i>E. coli</i> O157:H7: 560 CFU <i>E. coli</i> : 3,794 CFU	H1	17.19	35.26	Positive	17.20	35.38	Positive	17.30	36.06		Positive	
H2			17.11	35.52	Positive	17.33	34.88	Positive	17.39	34.10	Positive			
H3			17.12	35.01	Positive	17.36	35.39	Positive	17.20	34.31	Positive			
<i>E. coli</i> O157:H7: 6 CFU <i>E. coli</i> : 379 CFU		L1	18.67	34.01	Positive	18.75	34.18	Positive	18.70	33.70	Positive			
		L2	19.36	33.59	Positive	19.53	33.23	Positive	19.54	33.45	Positive			
		L3	19.36	34.01	Positive	19.25	33.65	Positive	19.42	34.01	Positive			

FAM (—), *E. coli* O157:H7 detection; HEX (—), internal control detection. BPW, buffered peptone water; CFU, colony forming units; STEC, Shiga toxin–producing *E. coli*.

Table 3. iQ-Check STEC VirX PCR Kit results for onions.

Matrix	Inoculum	Sample	Rep 1				Rep 2				Rep 3				PCR Curves		
			FAM	Cy5	HEX	Result	FAM	Cy5	HEX	Result	FAM	Cy5	HEX	Result			
Onions, fresh chopped 18 hr incubation BPW	No inoculum	0	N/A	N/A	32.15	Negative	N/A	N/A	31.94	Negative	N/A	N/A	32.44	Negative			
	<i>E. coli</i> O157:H7: 560 CFU <i>E. coli</i> : 3,794 CFU	H1	18.77	18.24	27.71	Positive	18.63	18.07	26.83	Positive	18.53	17.96	27.19	Positive			
		H2	18.72	18.18	28.25	Positive	18.54	18.03	27.01	Positive	18.39	17.69	26.86	Positive			
		H3	18.22	17.79	28.01	Positive	18.28	17.81	27.29	Positive	18.29	17.62	N/A	Positive			
	<i>E. coli</i> O157:H7: 6 CFU <i>E. coli</i> : 379 CFU	L1	20.63	19.9	30.26	Positive	20.53	19.9	29.97	Positive	20.99	20.05	29.64	Positive			
		L2	23.35	22.67	30.09	Positive	23.70	22.9	30.08	Positive	23.56	22.77	30.03	Positive			
		L3	20.94	20.42	30.82	Positive	20.73	20.29	30.67	Positive	20.63	20.23	30.33	Positive			
	Onions, fresh chopped 18 hr incubation BPW + STEC Supplement	No inoculum	0	N/A	N/A	32.19	Negative	N/A	N/A	32.06	Negative	N/A	N/A	31.86		Negative	
		<i>E. coli</i> O157:H7: 560 CFU <i>E. coli</i> : 3,794 CFU	H1	18.23	17.51	26.39	Positive	17.93	17.26	25.89	Positive	18.01	17.39	30.91		Positive	
H2			17.94	17.22	27.46	Positive	18.11	17.41	26.79	Positive	18.12	17.46	N/A	Positive			
H3			17.85	17.27	26.45	Positive	17.83	17.28	33.43	Positive	17.97	17.30	26.20	Positive			
<i>E. coli</i> O157:H7: 6 CFU <i>E. coli</i> : 379 CFU		L1	19.68	19.08	29.82	Positive	19.77	19.17	29.85	Positive	19.64	19.15	29.93	Positive			
		L2	20.37	19.52	30.14	Positive	20.42	19.73	30.18	Positive	20.14	19.55	29.49	Positive			
		L3	20.14	19.5	30.08	Positive	20.16	19.28	30.14	Positive	20.38	19.57	30.13	Positive			

FAM (—), *stx1/stx2* detection; Cy5 (—), *eae* detection; HEX (—) internal control detection. BPW, buffered peptone water; CFU, colony forming units; Cy, cyanine; STEC, Shiga toxin–producing *E. coli*.

Table 4. dd-Check STEC ddPCR results for onions.

Matrix	Inoculum	Sample	Accepted Droplets	stx Value, copies	eae Value, copies	Linkage, %	Sample Result	
Onions, fresh chopped 18 hr incubation BPW with DTT 1:100 dilution	No inoculum	0	19,679	0	0		Negative	
	<i>E. coli</i> O157:H7: 560 CFU	H1	17,108	16,406	16,014	88	Positive: linked targets	
		H2	18,459	35,281	32,900	77.8	Positive: linked targets	
		H3	18,982	35,492	32,780	75	Positive: linked targets	
	<i>E. coli</i> O157:H7: 6 CFU	L1	17,296	11,308	10,989	85.5	Positive: linked targets	
		L2	18,795	791	730	82.3	Positive: linked targets	
		L3	18,009	7,683	6,928	78.6	Positive: linked targets	
	Onions, fresh chopped 18 hr incubation BPW with DTT 1:1,000 dilution	No inoculum	0	19,549	0	0		Negative
		<i>E. coli</i> O157:H7: 560 CFU	H1	19,866	3,216	3,105	87.7	Positive: linked targets
H2			18,995	2,951	2,759	79.7	Positive: linked targets	
H3			17,869	3,592	3,176	75.4	Positive: linked targets	
<i>E. coli</i> O157:H7: 6 CFU		L1	18,786	1,146	1,084	86.5	Positive: linked targets	
		L2	18,042	64	66.6	74	Positive: linked targets	
		L3	16,675	709	689	78.1	Positive: linked targets	

BPW, buffered peptone water; CFU, colony forming units; DTT, dithiothreitol.

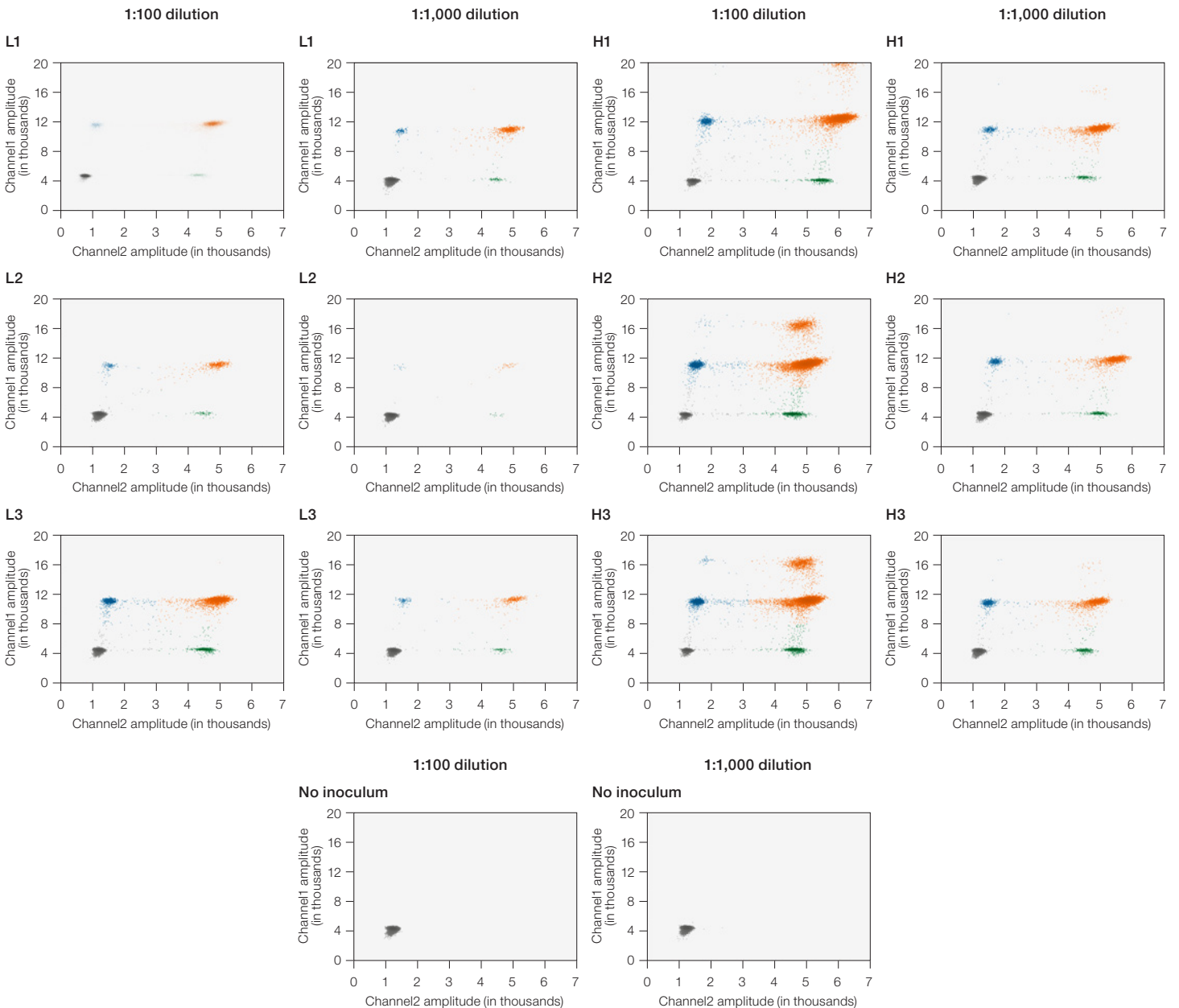


Fig. 1. 2-D amplitude graph of *E. coli* O157:H7 in onions when tested with dd-Check STEC Solution. The plots show droplets negative for both targets (■), droplets positive for *stx* only (■), droplets positive for *eae* only (■), and droplets positive for both *stx* and *eae* on the same genome (■).

Table 5. iQ-Check *E. coli* O157:H7 PCR Kit results for carrots.

Matrix	Inoculum	Sample	Rep 1			Rep 2			Rep 3			PCR Curves
			FAM	HEX	Result	FAM	HEX	Result	FAM	HEX	Result	
Carrots, fresh sliced 18 hr incubation BPW	No inoculum	0	N/A	33.20	Negative	N/A	33.08	Negative	N/A	32.52	Negative	
	<i>E. coli</i> O157:H7: 618 CFU	H1	17.83	N/A	Positive	17.88	38.94	Positive	17.82	38.35	Positive	
		H2	17.93	38.07	Positive	17.94	39.68	Positive	17.76	39.47	Positive	
		H3	18.53	37.54	Positive	18.27	38.40	Positive	18.04	38.19	Positive	
	<i>E. coli</i> O157:H7: 6 CFU	L1	20.08	36.82	Positive	19.95	37.48	Positive	19.78	35.84	Positive	
		L2	20.59	36.45	Positive	20.57	35.36	Positive	19.96	36.62	Positive	
		L3	19.93	36.49	Positive	20.02	N/A	Positive	19.34	39.71	Positive	

FAM (—), *E. coli* O157:H7 detection; HEX (—), internal control detection. BPW, buffered peptone water; CFU, colony forming units.

Table 6. iQ-Check STEC VirX PCR Kit results for carrots.

Matrix	Inoculum	Sample	Rep 1			Rep 2			Rep 3			PCR Curves			
			FAM	Cy5	HEX	Result	FAM	Cy5	HEX	Result	FAM		Cy5	HEX	Result
Carrots, fresh sliced 18 hr incubation BPW	No inoculum	0	N/A	N/A	32.04	Negative	N/A	N/A	32.11	Negative	N/A	N/A	31.85	Negative	
	<i>E. coli</i> O157:H7: 618 CFU	H1	19.3	18.05	31.4	Positive	19.14	18	31.16	Positive	19.12	18.02	31.20	Positive	
		H2	19.17	18.21	31.08	Positive	19.2	18.15	31.24	Positive	19.36	18.43	31.83	Positive	
		H3	19.51	18.65	30.76	Positive	19.41	18.51	31.10	Positive	19.36	18.32	32.85	Positive	
	<i>E. coli</i> O157:H7: 6 CFU	L1	20.94	20.24	31.23	Positive	20.77	20.12	30.82	Positive	20.48	19.87	31.75	Positive	
		L2	21.48	21.06	30.97	Positive	21.55	21.16	30.7	Positive	21.37	20.74	31.91	Positive	
		L3	21.1	20.27	30.56	Positive	21.17	20.35	31.08	Positive	20.91	19.80	31.80	Positive	

FAM (—), *stx1/stx2* detection; Cy5 (—), *eae* detection; HEX (—) internal control detection. BPW, buffered peptone water; CFU, colony forming units; Cy, cyanine.

Table 7. iQ-Check *Salmonella* PCR Kit results for cucumbers.

Matrix	Inoculum	Sample	Rep 1			Rep 2			Rep 3			PCR Curves
			FAM	HEX	Result	FAM	HEX	Result	FAM	HEX	Result	
Cucumbers, fresh sliced 18 hr incubation BPW	No inoculum	0	N/A	33.07	Negative	N/A	33.08	Negative	N/A	33.20	Negative	
	<i>S. Typhimurium</i> : 534 CFU	H1	21.04	34.4	Positive	21.15	33.62	Positive	21.01	33.29	Positive	
		H2	21.46	33.61	Positive	21.22	33.58	Positive	21.03	32.70	Positive	
		H3	21.36	33.49	Positive	21.18	33.17	Positive	21.15	32.98	Positive	
	<i>S. Typhimurium</i> : 7 CFU	L1	24.50	32.82	Positive	24.61	33.02	Positive	24.57	32.30	Positive	
		L2	23.82	33.06	Positive	23.68	32.55	Positive	23.46	32.75	Positive	
		L3	24.47	32.59	Positive	24.51	32.35	Positive	24.52	32.38	Positive	

FAM (—), *Salmonella* spp. detection; HEX (—) internal control detection. BPW, buffered peptone water; CFU, colony forming units.

Detection of *E. coli* O157:H7 in Carrots

When tested in triplicate with the iQ-Check *E. coli* O157:H7 PCR Kit, all inoculation levels showed positive results after 18 hr of incubation (Table 5). All inoculation levels also yielded positive results after 18 hr of incubation when tested with the iQ-Check STEC VirX PCR Kit (Table 6).

Detection of *Salmonella* in Cucumbers

When tested in triplicate with the iQ-Check *Salmonella* PCR Kit, positive results after 18 hr of incubation were obtained for all inoculation levels, with or without STEC Supplement (Table 7).

Conclusion

Onions, carrots, and cucumbers tested with iQ-Check and dd-Check PCR Kits were verified at the target enrichment ratio and incubation time for *E. coli* O157:H7 (onions and carrots) and *Salmonella* (cucumbers). Selective STEC Supplement was not required for the enrichment of onions. The dd-Check STEC Kit was able to determine that the strain used to inoculate onions contained both *stx* and *eae* virulence genes on the same genome. Samples were detected after 18 hr of incubation, which allows for a rapid response in the event of a contamination event.

Further Reading

- Center for Disease Control and Prevention. *E. coli* Outbreak Linked to Onions Served at McDonald's. [cdc.gov/ecoli/outbreaks/e-coli-O157.html](https://www.cdc.gov/ecoli/outbreaks/e-coli-O157.html) accessed January 13, 2025.
- Center for Disease Control and Prevention. *E. coli* Outbreak Linked to Organic Carrots. [cdc.gov/ecoli/outbreaks/e-coli-o121.html](https://www.cdc.gov/ecoli/outbreaks/e-coli-o121.html) accessed January 13, 2025.
- Center for Disease Control and Prevention. *Salmonella* Outbreak Linked to Cucumbers. [cdc.gov/salmonella/outbreaks/cucumbers-11-24/index.html](https://www.cdc.gov/salmonella/outbreaks/cucumbers-11-24/index.html) accessed January 13, 2025.
- Center for Disease Control and Prevention. Foodborne Illness Source Attribution Estimates – United States, 2022. [cdc.gov/ifsac/php/data-research/annual-report-2022.html](https://www.cdc.gov/ifsac/php/data-research/annual-report-2022.html) accessed January 13, 2025.

Visit [bio-rad.com/FoodSafety](https://www.bio-rad.com/FoodSafety) for more information.

BIO-RAD, DDP, DROPLET DIGITAL, DROPLET DIGITAL PCR, and QX200 are trademarks of Bio-Rad laboratories, Inc. in certain jurisdictions. IQ-CHECK is a trademark of Bio-Rad Europe GmbH in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2025 Bio-Rad Laboratories, Inc.

Purchase of Digital PCR and/or Single-Cell NGS Sample Preparation products (the "Products") from Bio-Rad Laboratories is subject to Bio-Rad Laboratories, Inc. Standard Terms and Conditions of Sale, which can be accessed at <https://www.bio-rad.com/en-us/terms-conditions>. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for you to distribute or resell the Products. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for the development or commercialization of diagnostic assays for use with the Products without a license from Bio-Rad. It is the user's obligation to obtain a commercial license from Bio-Rad for (i) all commercial uses (not just diagnostic uses) and (ii) sale of assays for use on Bio-Rad's dPCR and ddSEQ instruments. The Products and/or their use are covered by U.S. and foreign patents and/or pending patent applications owned by or under license to Bio-Rad Laboratories, Inc. See <https://www.bio-rad.com/en-us/trademarks>.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website [bio-rad.com](https://www.bio-rad.com) USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399 Canada 1 800 361 1808 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 080 007 7373 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23