

Advance Your Water Testing Capabilities: Molecular and Microbiological Approaches

Detect, Enumerate, and Identify Waterborne Pathogens and Quality Indicators

Common Threats in Water Systems

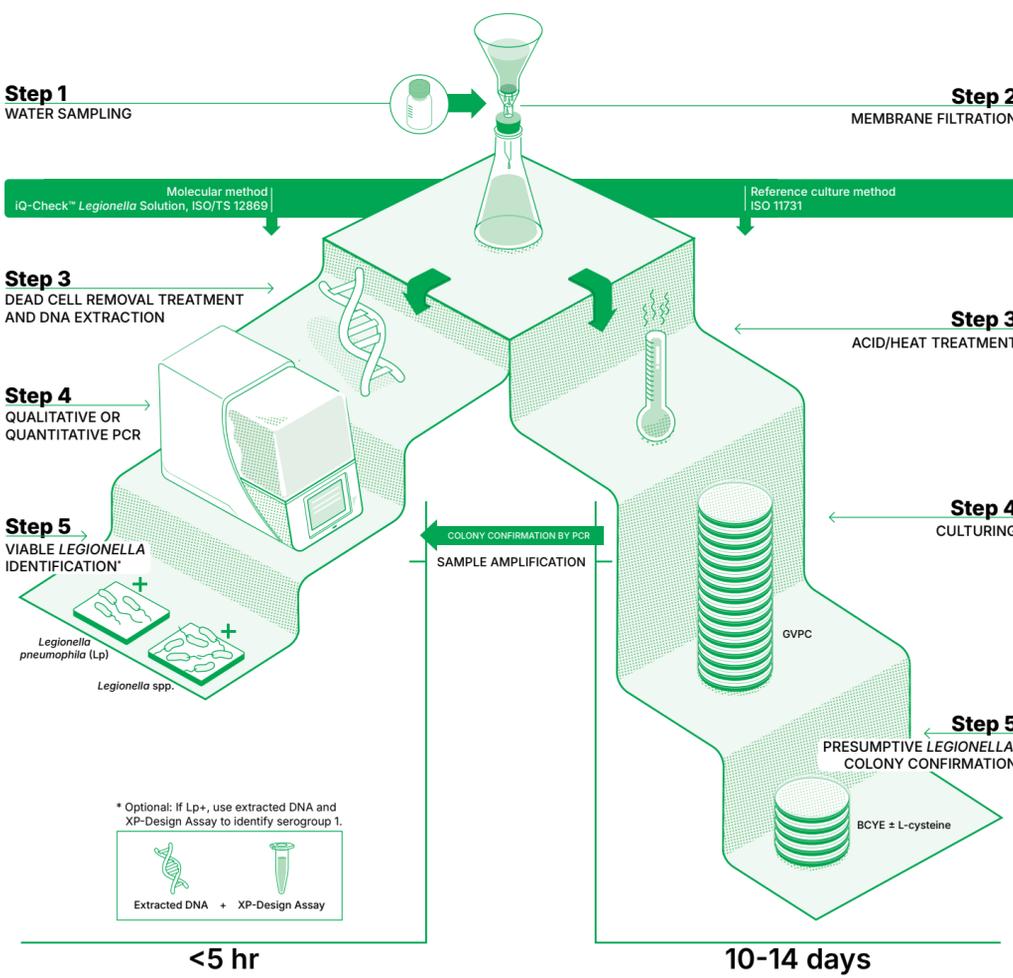
Pathogens and Indicators

Modern water testing combines molecular and culture-based methods to detect quality indicators and dangerous pathogens that pose significant health risks. These methods help identify the source, understand the risk, and guide targeted interventions.

Microorganism	Risk	Sources
<p><i>Legionella</i></p>	<ul style="list-style-type: none"> Causes Legionnaires' disease and Pontiac fever, especially in vulnerable groups Building owners and operators are responsible for prevention and control Thrives in biofilms within plumbing systems Spreads via inhaled water droplets 	
<p><i>E. coli</i> and coliforms</p>	<ul style="list-style-type: none"> Are indicators of potential fecal contamination Pose health risks to vulnerable populations Can enter systems through cracks, leaks, or runoff 	
<p><i>Pseudomonas aeruginosa</i></p>	<ul style="list-style-type: none"> Common in healthcare and industrial water systems Forms biofilms, making elimination difficult Thrives in moist environments like old plumbing, filters, and bottled water systems 	

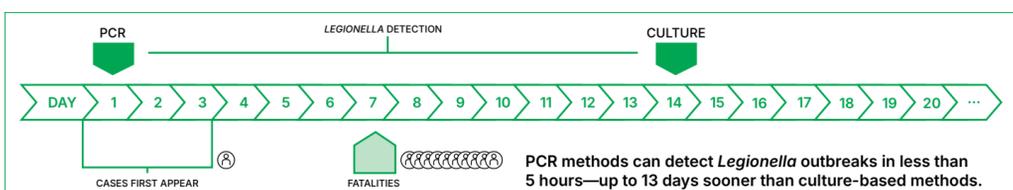
Detect *Legionella* Faster with PCR

Comparing Molecular and Microbiological Methods



Faster *Legionella* Testing Results Means Earlier Responses

PCR testing identifies contamination in less than 5 hr, while the culture method can take up to 14 days for results. In a *Legionella* outbreak, cases typically begin to appear around days 1–3, with fatalities occurring by day 7. Faster detection with PCR enables quicker response, improving crisis management and public health outcomes.



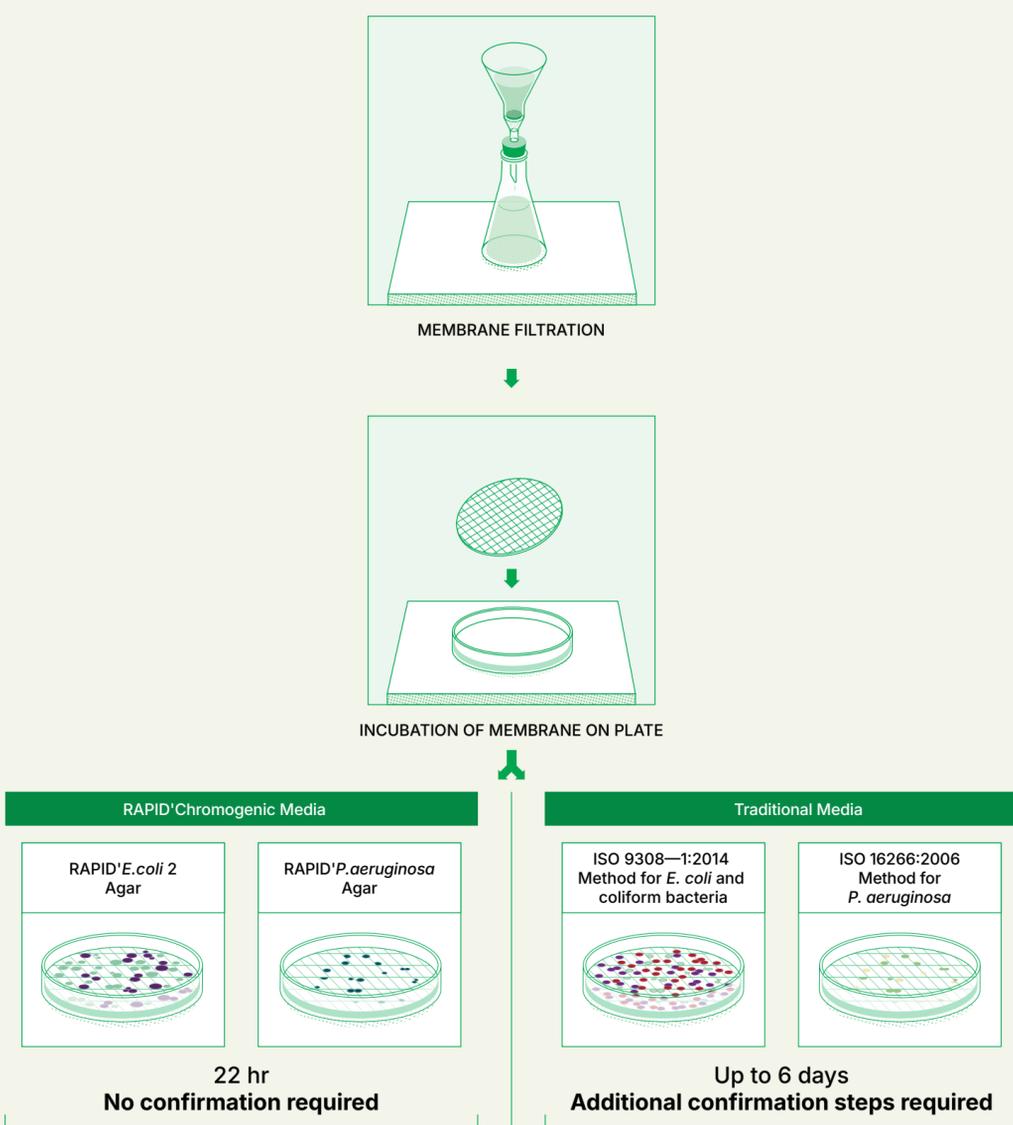
Stopping Outbreaks Before They Start: The Power of the iQ-Check *Legionella* Solution

Certified, reliable technology	Rapid, actionable results	Enhanced decision-making
AFNOR certified (ISO/TS 12869)	Same-day results	Qualitative and quantitative data
FDRS technology—detects only live cells	Quick screening of negative samples	Identifies Lp serogroup 1 with XP-Design Assay
Detects VBNC (viable but non-culturable) cells	Accelerate ISO 11731 workflows with PCR confirmation of presumptive colonies	Faster evaluation of remediation efforts

The iQ-Check *Legionella* Solution offers a rapid tool to complement culture methods.

Streamline Detection and Enumeration of Quality Indicators

Testing Methods for Quality Indicators



RAPID'Chromogenic Media provides results faster and with fewer steps.

Enhancing Quality Control with Chromogenic Media: Key Advantages for Water Testing

<p>RAPID'<i>E. coli</i> 2</p> <ul style="list-style-type: none"> US EPA and AFNOR validated alternative to ISO 9308 method Direct enumeration of <i>E. coli</i> and coliforms in one plate Results within 18 hr (AFNOR) or 22 hr (EPA) Easy interpretation: Excellent colony contrast coloration No confirmation step required 	<p>RAPID'<i>P.aeruginosa</i></p> <ul style="list-style-type: none"> AFNOR validated alternative to ISO 16266 method Direct enumeration of <i>Pseudomonas aeruginosa</i> Faster TTR: Results within 22 hr Easy interpretation: Excellent colony contrast coloration No confirmation step required
---	---

RAPID'Chromogenic Media offers a validated method for easy enumeration with faster results than traditional media.