



Reduce False Positives with Probe-Based Droplet Digital PCR *Mycoplasma* Testing



Testing for *Mycoplasma* contamination is a crucial step in the quality control process that requires sensitivity, specificity, and reproducibility. However, the standard SYBR® based quantitative PCR (qPCR) methods face challenges due to a lack of specificity and can result in high false-positive rates. The Bio-Rad VeriCheck Droplet Digital PCR (ddPCR) *Mycoplasma* Detection Kit provides the sensitivity you need via probe-based chemistry and offers the advantage of ddPCR inhibitor tolerance. Detect up to 112 *Mycoplasma* species and minimize false positives to avoid disruption of your development and manufacturing processes.

The VeriCheck ddPCR *Mycoplasma* Detection Kit provides:

- Highly specific results and low cross-reactivity with related species
- Reproducible results, reducing plate-to-plate and user-to-user variability
- Automated data analysis using positive control
- U.S. FDA 21 CFR Part 11 compatible QX Manager Software, Regulatory Edition (#12012172), offering audit trails with tracked protocol changes

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



BIO-RAD

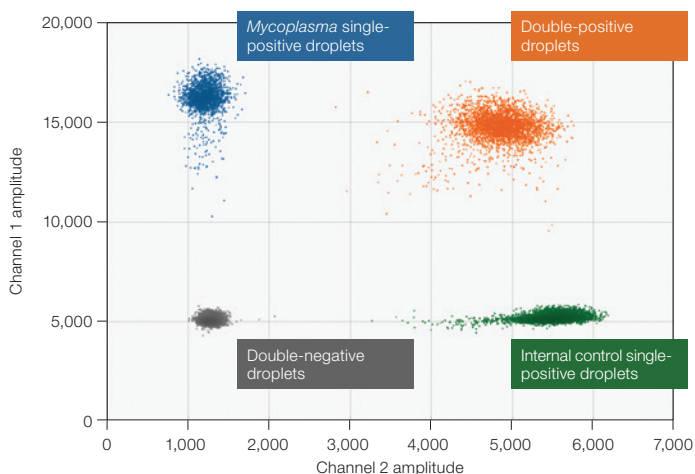
Limit of Detection (LOD)

Colony forming units (CFU)/ml	≤1
Genomic copies (GC)/reaction	≤6

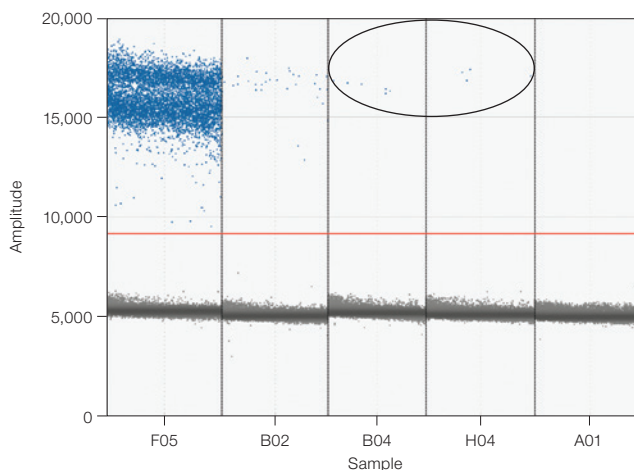
Nine *Mycoplasma* species at ≥95% of the replicate wells were positive for *Mycoplasma* extracted at 1 CFU/ml. The LOD in CFU/ml is determined to be 1 CFU/ml.

LOD in GC/well is determined to be ≤6 GC/well.

Performance Data



2-D amplitude plot for positive control well. Plot showing *Mycoplasma* signal in channel 1 (FAM) and internal control signal in channel 2 (HEX). Plot shows four droplet clusters. The gray cluster represents droplets that are double-negative for *Mycoplasma* and internal control, the blue cluster represents droplets that are positive for *Mycoplasma* only, the green cluster represents droplets that are positive for internal control only, and the orange cluster represents droplets that are double-positive for *Mycoplasma* and internal control.



Positive control based autothresholding. Positive control based autothresholding allows for accurate thresholding of an entire plate simultaneously. Accurate thresholds are applied to *Mycoplasma* samples with concentrations spanning the ddPCR dynamic range. Positive control used to threshold *Mycoplasma* samples (well F05). Samples at 35 genome copies (GC)/well (B02), 5 GC/well (B04), 3 GC/well (H04), and no template control (A01).

Ordering Information

Catalog #	Description
12013126	VeriCheck ddPCR <i>Mycoplasma</i> Detection Kit

BIO-RAD, DDP-PCR, and DROPLET DIGITAL PCR are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. SYBR is a trademark of Thermo Fisher Scientific Inc. All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website 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 1 905 364 3435 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 2 3473 4460 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

