

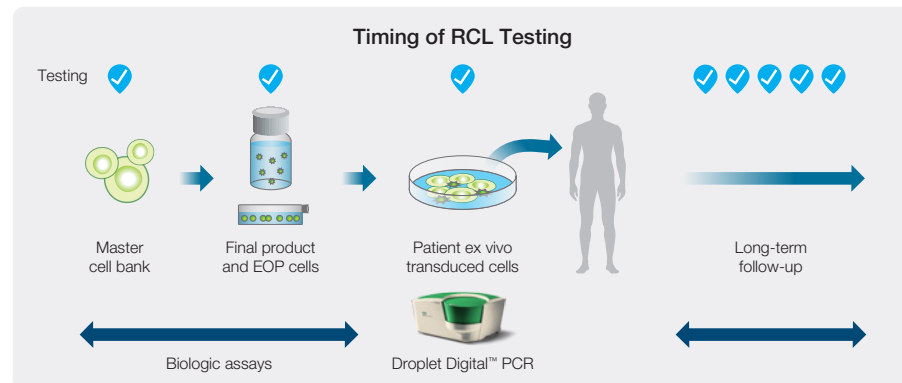
Vericheck ddPCR™ Replication Competent Lentivirus (RCL) Kit

A Faster, Quantitative, and Orthogonal Method for Testing RCL



Lentiviruses are commonly used as viral vectors in cell and gene therapy production, delivering and steadily integrating desired genes into the genome of the host cell. However, they may also present health hazards to humans due to integration-mediated transformation and the creation of a RCL that can infect nontarget cells. The U.S. Food and Drug Administration recommends testing sufficient supernatant to ensure a 95% probability of RCL detection if present at a concentration of more than 1 RCL/dose equivalent. While the current regulatory guidelines recommend a cell culture–based method for RCL testing, it is cumbersome and takes 30–45 days from sample to results.

Bio-Rad's Vericheck ddPCR RCL Kit is fast, cost-effective, specific, and a fully validated solution for lentivirus testing.



EOP, end of production.

Quick turnaround time



Get results in <8 hours, compared to 30–45 days for culture-based methods

Qualitative and quantitative results



Perform automated data analysis with the ability to use prevalidated assay protocol files (APF)

Substantial savings



Economical alternative to culture-based methods

Specific and sensitive

99.9% Specificity

Limit of detection: 0.87 copies/μl
≥95% probability of detection

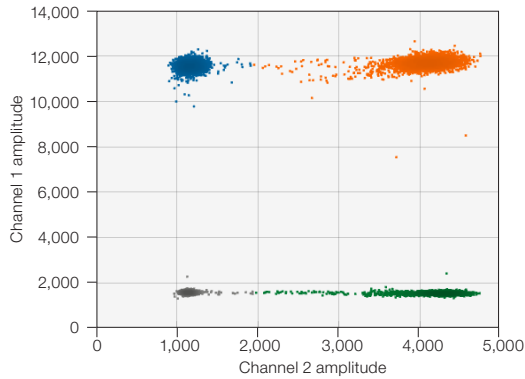


Fig. 1. 2-D amplitude plot for the VeriCheck ddPCR Replication Competent Lentivirus (RCL) Kit. The plot shows the RCL signal in channel 1 (FAM) and the internal control signal in channel 2 (HEX). The blue cluster is single-positive for RCL DNA, the green cluster is single-positive for internal control DNA, the orange cluster is double-positive for both RCL and internal control DNA, and the gray cluster is double-negative.

The ddPCR RCL Kit Includes:

- ddPCR Supermix for Residual DNA Quantification, 2x supermix
- ddPCR RCL Internal Control
- ddPCR RCL Positive Control
- Nuclease-Free Water (Negative Control)
- ddPCR RCL Assay

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

BIO-RAD, DDPCR, DROPLET DIGITAL, and DROPLET DIGITAL PCR are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2023 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** 1 905 364 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** 82 080 007 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

