

## ZDC Multiplex RT-PCR Assay

Bio-Rad Laboratories introduces the ZDC Multiplex RT-PCR Assay for the simultaneous detection of Zika (ZIKV), dengue (DENV 1–4), and chikungunya (CHIKV) viral RNAs by real-time PCR. Benefiting from a robust one-step RT-PCR protocol and expertly designed primers and probes, this kit enables sensitive identification of viral targets together with multiplex detection of internal and sample controls.

- Wet-lab validated on the CFX96 Touch™ Real-Time PCR Detection System and Applied Biosystems' 7500 Real-Time PCR System and compatible with most qPCR systems on the market
- Compatible with RNA extracted from multiple starting materials (serum, plasma, serum-free body fluids, cerebrospinal fluids, and mosquitoes)
- High-standard GMP manufacturing procedures used for key components of the kit
- All controls required to validate the results included in the kit
- Compatible with RNase P Detection Assays for further expansion of the range of controls

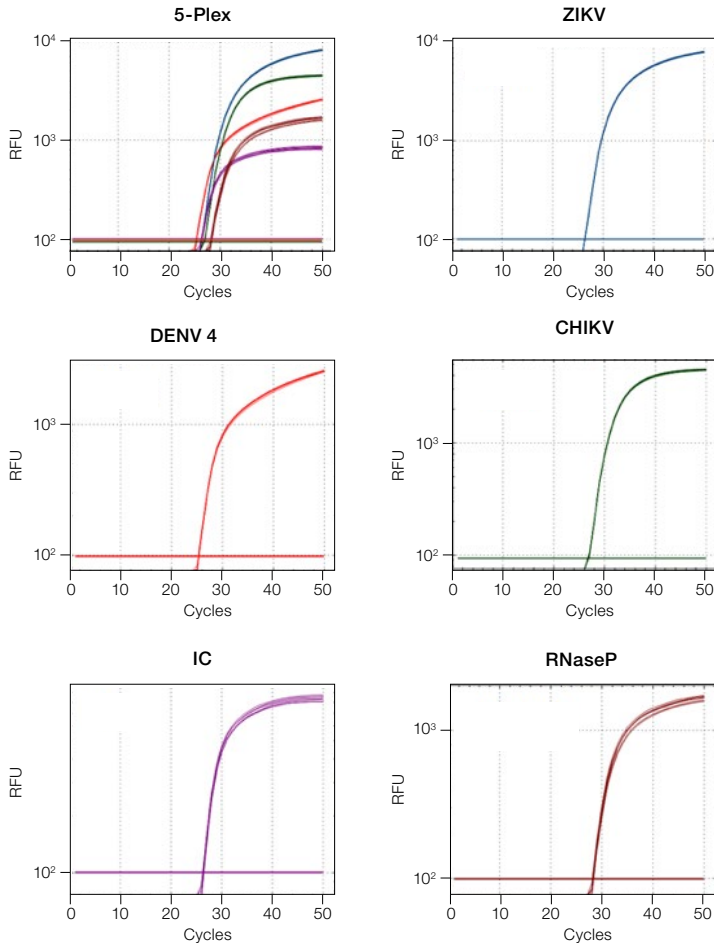
Visit [bio-rad.com/ZDC](https://www.bio-rad.com/ZDC) for more information.



## ZDC Multiplex RT-PCR Assay

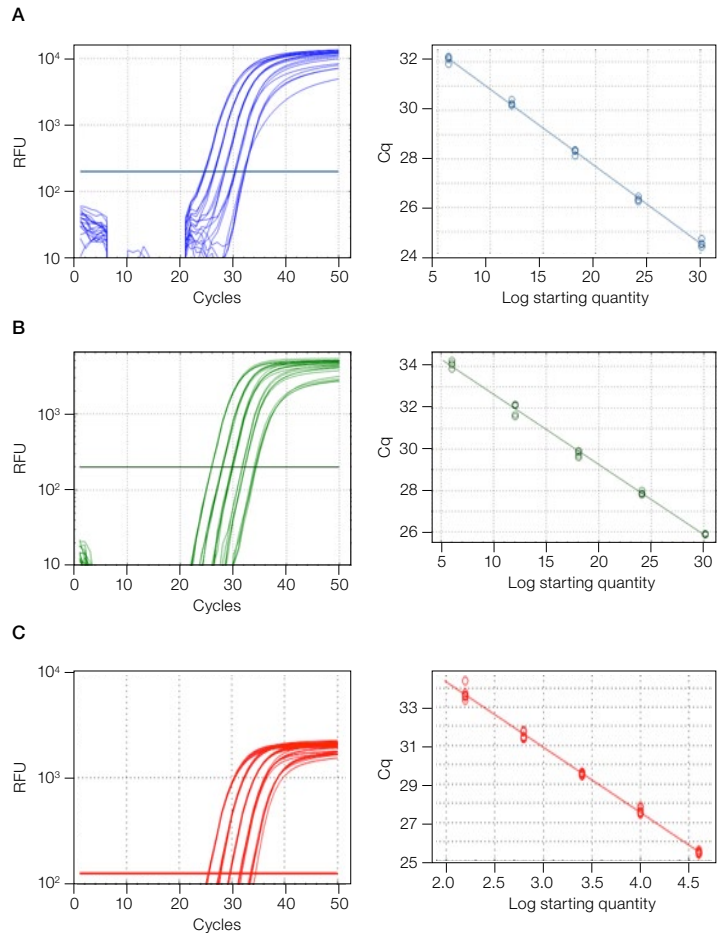
Bio-Rad's ZDC Multiplex RT-PCR Assay is a wet-lab validated, RT-PCR multiplex assay for the detection of Zika, dengue, and chikungunya viral RNA from human serum, mosquito tissues, and more. Including GMP-manufactured key components and all the controls required to validate your results, the ZDC Assay provides researchers with a sensitive and robust tool for the simultaneous detection of these arboviruses.

### One-step 5-plex capability



**One-step 5-plex capability.** ZDC Multiplex RT-PCR Assay, including RT-PCR reactions for the five potential targets: ZIKV, DENV 4, CHIKV, internal control (IC), and RNase P. The experiment was performed on the CFX96 Touch Real-Time PCR System.

### Excellent linearity and efficiency



**Excellent linearity (>99% R<sup>2</sup>) and efficiency (>98%) for all targets.** Five-point (1:4) serial dilution on the CFX96 Touch Real-Time PCR System of ZIKV (A), DENV 2 (B), and CHIKV (C).

### Ordering Information

Catalog #	Description
12003818	<b>ZDC Multiplex RT-PCR Assay</b> , real-time PCR assay for the detection of Zika, dengue, and chikungunya viral RNAs (200 reactions)
12004601	<b>RNase P Assay (Cy5.5)</b> , pkg of 1, vial, 440 µl of 12.5x concentrated primers/probe solution
12004602	<b>RNase P Assay (TAMRA)</b> , pkg of 1, vial, 440 µl of 12.5x concentrated primers/probe solution
12004860	<b>ZDC Clear Control Solution</b> , pkg of 1, 5 ml tube, for use with RNase P Assay and ZDC Multiplex RT-PCR Assay
12005062	<b>ZDC Buffer T</b> , pkg of 5, 1.8 ml tubes, for use with ZDC Multiplex RT-PCR Assay

The ZDC Multiplex RT-PCR Assay is intended for research use only and not for use in diagnostic procedures.

Applied Biosystems is a trademark of Thermo Fisher Scientific. Cy is a trademark of GE Healthcare. TAMRA is a trademark of Life Technologies Corporation.

Bio-Rad's thermal cyclers and real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00 **France** 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026 674 55 05 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

