



## ddPCR Assays for Adeno-Associated Virus (AAV) Vector Genome Titer



Bio-Rad Droplet Digital PCR (ddPCR) Assays for AAV vector genome titer utilize the precision and sensitivity of ddPCR technology to deliver an absolute count of AAV vector genome copies in your sample. These assays work seamlessly with the Bio-Rad ddPCR product portfolio to offer the convenience of a complete workflow solution required to help you validate the quality of your therapeutic product with confidence.

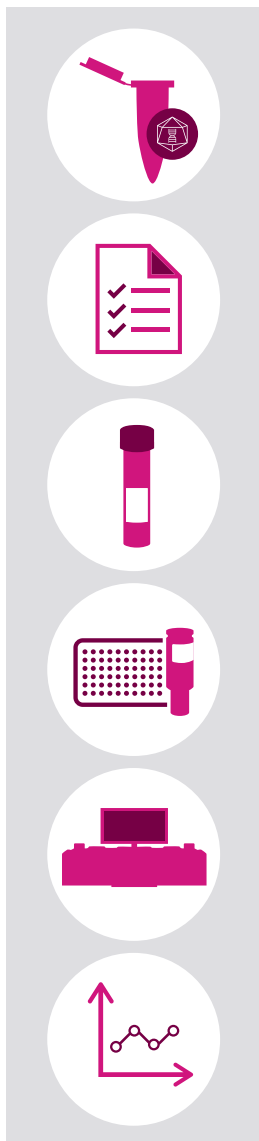
- Custom ddPCR Cell and Gene Therapy Assays generated by the Cell and Gene Therapy Design Engine
- Ready-to-order ddPCR Expert Design Assays for common vector and plasmid elements

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



**BIO-RAD**

## Complete ddPCR Workflow Solution for AAV Vector Genome Titer



### Assays

- ddPCR Cell and Gene Therapy Assays: custom assays for your gene of interest, available in 6 fluorophore choices
- Vector Backbone Assays: ddPCR Expert Design Assays for common viral vector backbone and plasmid elements

### Protocol

- Optimized protocol designed as a starting point for the most precise measurement of AAV vector genome titer
- Droplet Digital PCR offers a significantly lower coefficient of variation (CV <10%) compared to quantitative PCR across both in-process and purified samples

### Supermixes

- ddPCR Supermix for Probes (No dUTP): ready-to-use 2x mix for probe-based Droplet Digital PCR
- ddPCR Multiplex Supermix: ready-to-use 4x mix for multiplex probe-based Droplet Digital PCR on the QX ONE Droplet Digital PCR System

### Consumables

- Specially formulated Droplet Generation and Droplet Reader Oils unique to each ddPCR System
- Cartridges, 96-well plates, and other consumables unique to each ddPCR System

### Platforms

- QX ONE ddPCR System: fully integrated ddPCR workflow and 5-plate capacity, allowing for walk-away operation; 4 color channels for advanced multiplexing
- QX200 AutoDG Droplet Digital PCR System: automated droplet generation, eliminating user-to-user variability by providing more consistent droplet counts across users and plates

### Software

- QX ONE Software, Regulatory Edition: robust software package that works with the QX ONE ddPCR System and permits operation in compliance with U.S. FDA 21 CFR Part 11 regulations
- QX Manager Software, Regulatory Edition: works with the QX200 and QX200 AutoDG ddPCR Systems and permits operation in compliance with U.S. FDA 21 CFR Part 11 regulations

### Ordering Information

Speak with your Bio-Rad Sales Specialist to understand what a complete ddPCR AAV vector genome titer solution would look like for you.

BIO-RAD, DDPCR, DROPLET DIGITAL PCR, and QX ONE are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

Purchase of Digital PCR products from Bio-Rad Laboratories includes a limited, non-transferable right under our intellectual property for you to use the products for your internal research purposes only. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for you to distribute or resell the products or to use the products for commercial applications, such as contract services or fee for services. No rights are granted for diagnostic uses. Information concerning a license for such uses can be obtained from Bio-Rad Laboratories. It is the responsibility of the purchaser/end user to acquire any additional intellectual property rights that may be required. The QX200, QX200 AutoDG, and QX ONE Droplet Digital PCR Systems and/or their use is covered by claims of U.S. patents, and/or pending U.S. and non-U.S. patent applications owned by or under license to Bio-Rad Laboratories, Inc. See [bio-rad.com/en-us/trademarks](http://bio-rad.com/en-us/trademarks) for details.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Website** [bio-rad.com](http://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

