

## ddPCR™ Droplet Reader Oil EcoTank for Probes

High-Capacity Droplet Reader Oil Reservoir for Use with the QX200™ Droplet Reader



The Bio-Rad ddPCR Droplet Reader Oil EcoTank for Probes enables droplet reader oil recirculation for a lower cost and greener workflow while delivering the same best-in-class performance on the QX200 Droplet Digital™ PCR System.

### Features and Benefits



Run up to 50 plates with a single EcoTank



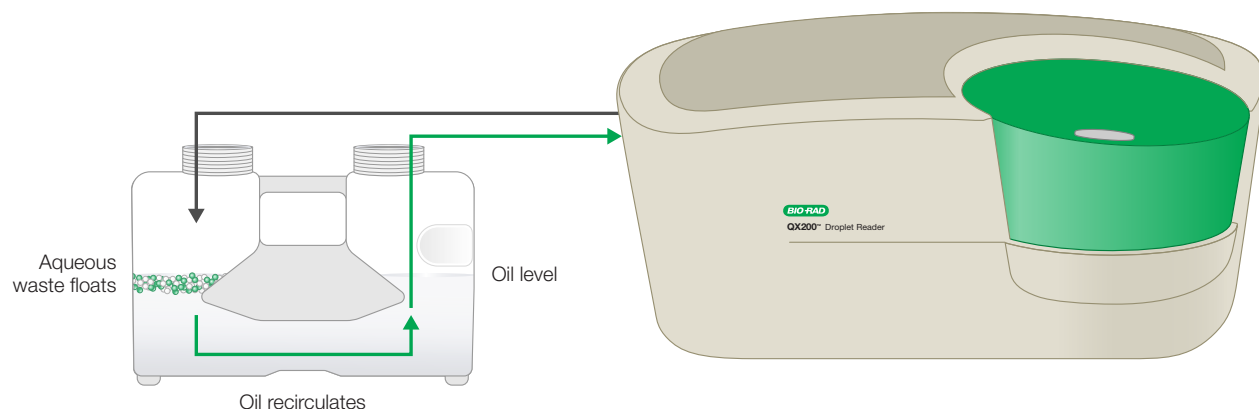
Reduce waste and per-reaction running cost



Implement the EcoTank with a simple software update



Achieve the same high-quality data you expect from the QX200 ddPCR System



**Fig. 1. Oil recirculation with the EcoTank.** The EcoTank separates floating aqueous waste generated by the ddPCR instrument, allowing droplet reader oil to be recirculated for up to 50 runs.

### How It Works

The ddPCR Droplet Reader Oil EcoTank for Probes is built for use during the droplet analysis step of the digital PCR workflow on the QX200 Droplet Reader and fits in the existing oil location without any hardware modification. The EcoTank distributes droplet reader oil and collects aqueous waste generated by the instrument into a single partitioned container (Figure 1). It ensures floating waste droplets remain on the waste side of the instrument for another droplet reading cycle. The EcoTank has been optimized and tested for probe-based detection.

	Droplet Reader Oil	
	Standard Bottle	EcoTank
Number of plates	8	50
Cost per well	\$\$\$	\$\$
Bottle change frequency*	Once every 2 weeks	Once every 10 weeks

\* Based on five plates per week.

### Ordering Information

Catalog #	Description
12019215	ddPCR Droplet Reader Oil EcoTank for Probes, 750 ml

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

BIO-RAD, DDPCR, DROPLET DIGITAL, and QX200 are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. Copyright © 2023 by 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**

**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

