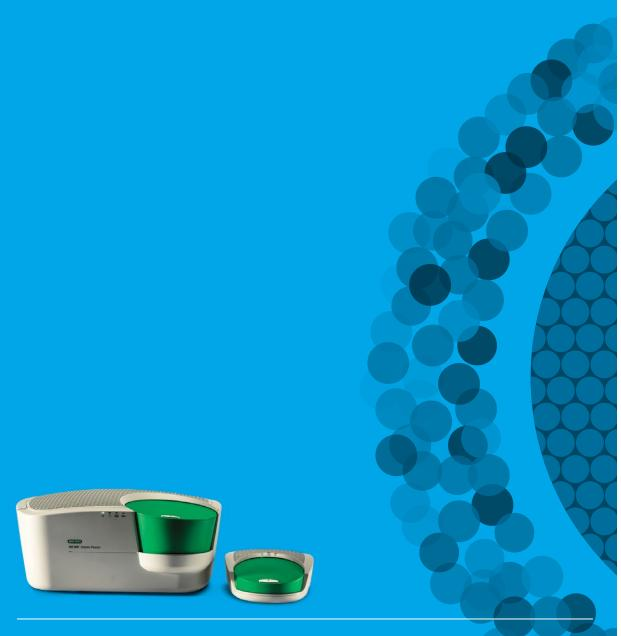


QX200[™] Droplet Digital[™] PCR System







Droplet Digital PCR (ddPCR[™]) is Bio-Rad's unique digital PCR technology. With unrivaled precision, Droplet Digital PCR provides absolute quantification of target DNA or RNA molecules without the use of standard curves. It addresses the lack of scalable and practical technologies for digital PCR implementation. The new QX200 Droplet Digital PCR System puts this powerful technology in your hands, ready to unveil new worlds of research at previously unattainable levels.

PCR becomes "digital" through sample partitioning and the subsequent statistical analysis of target detection across the partitions. The QX200 System uses advanced microfluidics technology to achieve partitioning on a massive scale, generating 20,000 highly uniform nanoliter-sized droplets per sample. For even higher sensitivity, it is possible to combine wells and analyze millions of droplets for a single sample.

Simple and reliable, Droplet Digital PCR technology has already led to several research breakthroughs in fields such as cancer biomarker discovery, infectious diseases, genomic alterations, and gene expression. This research is only the beginning. What discoveries will you make using Droplet Digital PCR?

A VERSATILE AND SCALABLE WORKFLOW



Droplet Digital PCR has a simple, user-friendly experiment setup designed to process eight samples at a time. The process easily scales up to run a 96-sample experiment in 5 hours with minimal hands-on time. When higher throughput is required, multiple 96-sample experiments can be run in a day.

Bio-Rad's QX200 ddPCR System combines water-oil emulsion droplet technology with microfluidics. The QX200 System consists of two instruments — a droplet generator and a droplet reader — and associated consumables. The droplet generator partitions each sample into 20,000 uniform nanoliter-sized droplets in which nucleic acid molecules are distributed in a random fashion. Each droplet serves to partition the reactions. Droplets are transferred to a 96-well PCR plate and PCR is performed to end point in a thermal cycler. The droplets stream single file through the reader for fluorescence analysis. Positive droplets, which contain at least one copy of the target DNA or RNA molecule, exhibit increased fluorescence compared to negative droplets. The fraction of PCR-positive droplets enables the target to be quantified according to the Poisson distribution.

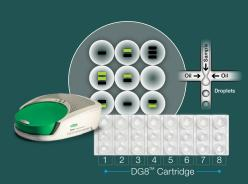
The QX200 System works with both hydrolysis probes and EvaGreen fluorescence detection chemistries, and its flexible design allows for high throughput and ultra-sensitive detection.



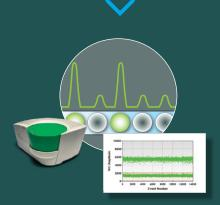
KEY BENEFITS

- Achieve absolute quantification without the use of a standard curve
- Design scalable assays for high sensitivity or high throughput
- Expand applications using flexible ddPCR chemistry EvaGreen or probes









Prepare ddPCR reaction mix

- Combine DNA sample, primers, and/or probes with one of Bio-Rad's ddPCR Supermixes
- Fully validated PrimePCR ddPCR Assays can be used

Generate droplets

- Load the ddPCR reaction mix into the wells of a droplet generator cartridge
- 8 x 20,000 droplets are generated from each run in the QX200 Droplet Generator
- Target DNA (—) and background DNA (—) are randomly distributed in droplets

Perform PCR with EvaGreen or hydrolysis probes

- Transfer the droplets to a 96-well PCR plate and seal the plate
- Run the PCR protocol

Read and analyze results

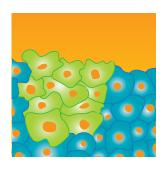
- After PCR, load the 96-well PCR plate into the QX200 Droplet Reader
- Positive and negative droplets in each sample are read
- Analyze concentrations with QuantaSoft[™] Software

QUANTIFYING BIOMARKERS FOR CANCER



The level of sensitivity offered by the QX200 ddPCR System in quantifying cancer biomarkers overcomes the limitations posed by other methods. Using Droplet Digital PCR technology, researchers are now able to observe finer quantitative distinctions among mutations, better identifying their potential roles in cancer.

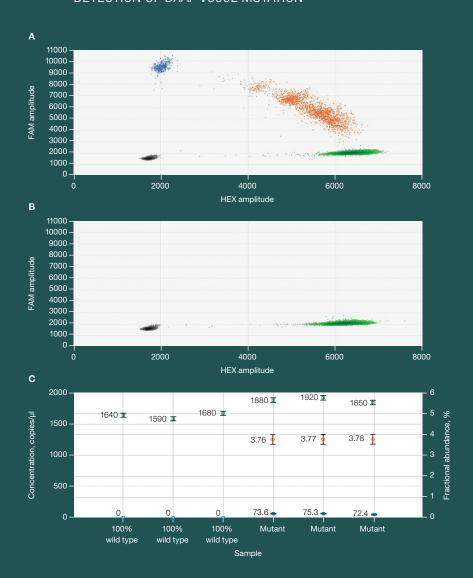
Cancer-associated mutations often evade detection due to their low concentrations relative to the background of wild-type DNA in a given sample. With its high sensitivity, the QX200 System can easily scale to quantify target concentrations as low as one out of 1,000,000 (0.0001%) total copies. What was previously undetectable with other methods can now be quantified with Droplet Digital PCR.



SUCCESS STORY

Detecting *T790M* mutation in epidermal growth factor receptor *(EGFR)*, an important therapeutic target in some lung cancers, allows for a better understanding of resistance to tyrosine kinase inhibitor therapies. However, other techniques lacked the sensitivity to reliably identify the *T790M* allele in a high background of *EGFR* wild type. Using Droplet Digital PCR, a researcher in the biotechnology industry was able to develop a highly accurate assay to quantify *T790M* concentration.

DETECTION OF BRAF V600E MUTATION



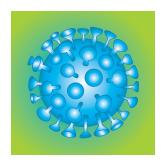
High sensitivity of the QX200 System allows quantification of BRAF V600E mutation in the presence of wild-type DNA using PrimePCR ddPCR Assays. A, 2-D fluorescence amplitude plot generated by QuantaSoft Software shows triplicate wells of a mixed mutant and wild-type sample. The black cluster on the plot represents the negative droplets, the green cluster represents the droplets that are positive for wild-type DNA only, the blue cluster represents the droplets that are positive for mutant DNA only, and the orange cluster represents the droplets that are positive for both mutant and wild-type DNA. B, 2-D fluorescence amplitude plot shows three replicates of a wild-type—only sample. C, fractional abundance plot shows the percentage frequency (orange markers) of the mutant DNA in a wild-type DNA background. The blue markers indicate the concentration of mutant DNA (copies/µl) and the green markers indicate the concentration of wild-type DNA (copies/µl) in each of three replicate samples. All error bars are generated by QuantaSoft Software and represent a

INSIGHT INTO VIRAL RESERVOIRS



Precisely quantifying viral load is crucial for characterizing disease states and developing and validating therapies. Viral reservoirs fluctuate, sometimes dropping to very low levels that are nonetheless significant. Obtaining exact and reproducible measurements of viral DNA or RNA often represents the difference between success and failure in studying many different pathogens.

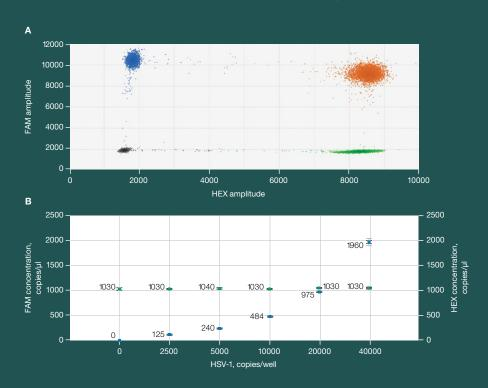
The QX200 System has the sensitivity to register extremely minute quantities of viral genetic material and distinguish it from complex mixtures. Investigators can obtain actual copy numbers of viral genome by generating complementary DNA through reverse transcription. Along with precision and reliability, the QX200 System also offers users the throughput needed to process samples with great efficiency.



SUCCESS STORY

A researcher compared quantitative PCR (qPCR) and Droplet Digital PCR approaches for detecting residual human immunodeficiency virus (HIV) infection in a clinical sample set, which included an infant who was functionally cured of HIV. The researcher found that Droplet Digital PCR was five times more sensitive than qPCR when measuring HIV DNA copies/million cells, and more than 20 times more accurate when measuring viral long terminal repeats (2-LTR circles).

DETECTION OF HERPES SIMPLEX VIRUS AND β2 MICROGLOBULIN



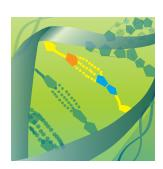
Droplet Digital PCR enables precise and reproducible detection of herpes simplex virus 1 (HSV-1) and 62 microglobulin (B2M) targets in a duplex assay. A, 2-D fluorescence amplitude plot shows three merged replicate samples of 10,000 copies/well of HSV-1 duplexed with 20,000 copies/well of B2M. The black cluster on the plot represents the negative droplets, the green cluster represents the droplets that are positive for HSV-1 only, the blue cluster represents the droplets that are positive for B2M only, and the orange cluster represents the droplets that are positive for both HSV-1 and B2M targets. B, concentration plot shows merged triplicate wells across an HSV-1 sample dilution series from 0 to 40,000 copies in a constant background of 20,000 copies (66 ng human genomic DNA) of B2M. The blue markers indicate HSV-1 copies/µl and the green markers indicate B2M copies/µl. All error bars are generated by QuantaSoft Software and represent a 95% confidence interval.

DISTINGUISHING GENOMIC VARIATIONS



Copy number variation (CNV) is a prominent source of interindividual variability in the human genome. CNV has also been associated with cancers, neurological and autoimmune diseases, and adverse drug responses. Reliably identifying such CNVs represents an important capability in cutting-edge research.

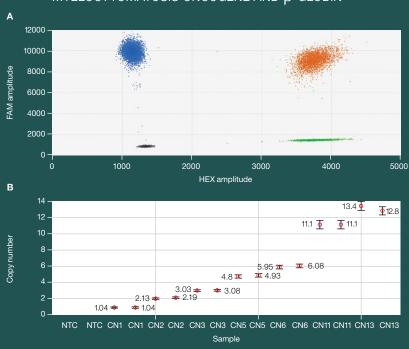
The major technical hurdle in copy number (CN) assessment is the ability to discriminate, with statistical confidence, between consecutive CN states. Fundamentally, as CN state increases, the percentage difference in target genomic material between states decreases, making it harder to measure CNV at higher orders. By partitioning the amplification reaction across thousands of droplets, the QX200 System can resolve high consecutive copy number states (such as five vs. six) with an excellent degree of statistical confidence.



SUCCESS STORY

A researcher investigating a highly duplicated gene involved in human brain size evolution needed a method to validate small CN variations in test samples. Droplet Digital PCR proved much more effective than other methods the researcher had previously used, showing greater precision in detecting small differences at high CN states, as well as excellent reproducibility. These improvements in CNV quantification will allow the researcher to ask more complex questions regarding neurogenetic disorders such as micro/macrocephaly, schizophrenia, and autism.

DETECTION OF COPY NUMBER FOR MYELOCYTOMATOSIS ONCOGENE AND β -GLOBIN



The QX200 System provides superior resolution with copy number (CN) calls from CN1 to CN13. A, 2-D fluorescence amplitude plot shows duplicate wells of a CN6 sample duplexed with myelocytomatosis oncogene (MYC) and β -globin (HBB) assays. The black cluster on the plot represents the negative droplets, the green cluster represents the droplets that are positive for MYC only, the blue cluster represents the droplets that are positive for HBB only, and the orange cluster represents the droplets that positive for both MYC and HBB targets. B, copy number plot shows samples ranging from CN1 to CN13 with precise duplicate well values. All error bars are generated by QuantaSoft Software and represent a 95% confidence interval. NTC, no template control.





Bio-Rad's ddPCR platform can be easily incorporated into the next-generation sequencing (NGS) library preparation workflow to precisely quantify and balance sequencing libraries on sequencers. In addition to precise quantification, the data plots generated by the QX200 System are rich with qualitative library information, which is not available with other currently available methodologies. Digital PCR enables consistent library loading and efficient utilization of next-generation sequencers. After the NGS run, Droplet Digital PCR can also be used to validate the results obtained by NGS, including genomic alterations such as single nucleotide polymorphisms, mutations, and copy number variations.





With its high sensitivity, the QX200 System is especially well-suited for discerning and quantifying rare transcripts, offering a better understanding of actual RNA function. With its high precision, the QX200 System can be used to detect small fold changes in expression of a target gene between samples, providing insights into the gene's physiological implications. The ability to measure absolute values of RNA molecules adds a new level of insight to gene expression studies. The QX200 ddPCR System allows for such measurements without requiring a standard curve.



SUCCESS STORY

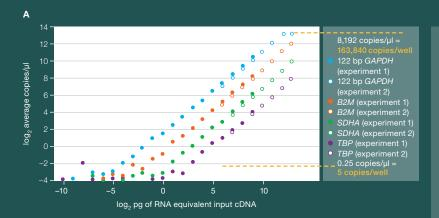
A research group used Droplet Digital PCR to accurately quantify five NGS libraries of varying concentration. Bio-Rad's ddPCR platform allowed the researchers to achieve an optimal cluster density of approximately 800,000/mm² required for sequencing on the Illumina MiSeq platform. Tests for library balancing showed that libraries could be balanced to within less than 15% of each other with a confidence interval of 95%.



SUCCESS STORY

A researcher used Droplet Digital PCR to study the differential expression of huntingtin (HTT) mRNA encoded by both wild-type and mutant HTT genes. Quantifying these different transcripts of RNA gave the researcher more insight into their cellular functions, and laid the groundwork for better understanding disease mechanisms.

VALIDATION FOR NEXT-GENERATION SEQUENCING



Droplet Digital PCR provides nearly 1,000-fold greater sensitivity and an amplification bias-free approach compared to RNA-Seq for validation. A, concentration plot shows a twofold dilution of cDNA generated using Bio-Rad's iScript™ Advanced cDNA Synthesis Kit for RT-qPCR. Two independent measurement sets were made: one at a high concentration range and one at a lower range, with four points overlapping. B, table shows comparative results between Droplet Digital PCR and MiSeq. Droplet Digital PCR detected thousands of copies per well with low-abundance transcripts (*TBP* and *GUSB* genes) using an input of 100 ng total RNA, which demonstrates enhanced sensitivity. The MiSeq Sequencer detected only single-digit reads per kilobase per

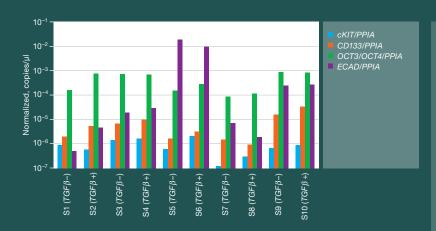
million reads (RPKM) for these transcripts.

В

Gene	Droplet Digital PCR, copies/well		MiSeq, RPKM			
	100 ng RNA 4 replicates	1,000 ng 4 replicates	100 ng 2 replicates		1,000 ng 2 replicates	
GAPDH	1,671 ± 115*	16,275 ± 479*	974	953	1,077	1,061
B2M	504 ± 46*	3,450 ± 155*	233	251	229	237
SDHA	139 ± 14*	1,131 ± 81*	51	50	59	65
HPRT1	15,781 ± 2,310	140,705 ± 11,059	34	22	25	25
TBP	3,650 ± 178	31,625 ± 1,010	3	3	6	3
GUSB	1,794 ± 53	15,731 ± 1,134	9	12	13	12

^{*} Obtained with cDNA diluted 200-fold because of its high abundance.

GENE EXPRESSION ANALYSIS OF STEM CELL MARKERS



High sensitivity and precision of the QX200 System allow reliable detection of small fold changes with rare transcripts. The $TGF\beta$ signaling pathway plays a role in ovarian stem cell regulation, and the addition of $TGF\beta$ will increase the number of stem cells in the population, increase the expression of putative stem cell markers, and decrease terminal differentiation markers like E-cadherin (ECAD). Data show expression levels of cKIT, CD133, OCT3/OCT4, and ECAD genes across ten samples with and without $TGF\beta$ treatment, normalized to PPIA in this study. Reproducible detection of cDNA copies as low as 0.5-1 target/ μ l in a $20~\mu$ l reaction was achieved. S, sample.

ABSOLUTE QUANTIFICATION WITH EvaGreen



Through droplet partitioning and statistical analysis, Droplet Digital PCR provides absolute quantification without the need for a standard curve, bringing substantial improvements to many applications that require detection of low-abundance target copies or distinction of small fold changes between samples. Combining Droplet Digital PCR with the power of EvaGreen detection chemistry enables genotyping and gene expression analysis for low-level target genes. The QX200 ddPCR System enables absolute quantification using EvaGreen chemistry, making such assays ideal for target DNA measurement, viral load analysis, and microbial quantification.

VALIDATED ddPCR ASSAYS



Bio-Rad's PrimePCR ddPCR* Assays are fully wet-lab validated assays, expertly designed to provide single-copy PCR resolution without the use of a standard curve. These mutation detection and copy number assays have been validated specifically for digital PCR.

Bio-Rad's PrimePCR Real-Time PCR Assays* can also be used on the QX200 Droplet Digital PCR System with QX200 ddPCR EvaGreen Supermix. Choose from more than 40,000 existing human and mouse gene expression assays.



SUCCESS STORY

Precision cancer medicine involves rapid assessment of cancer mutations, with certain genetic errors serving as indicators for a specific therapy. A researcher was challenged with the composition and degradation of clinical samples that were used for assessment of FLT3 mutations for gastric tumors (sporadic diffuse). By partitioning the samples' DNA templates, Droplet Digital PCR removed the effects of PCR inhibitors and bias in the data, resulting in a more reliable quantification of the mutations present in the samples.

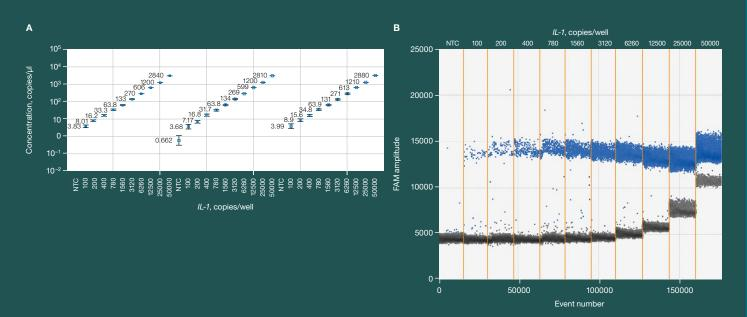


SUCCESS STORY

Designing an optimal probe assay may pose challenges and delay generation of reliable data. A researcher spent a great deal of time trying to design probes and primers for *EGFR L858R*, a receptor for epidermal growth factor, which is associated with lung tumors. Using the fully validated PrimePCR ddPCR Mutation Detection Probe Assay for *EGFR L858R* eliminated these tedious optimization steps, allowing the researcher to focus on the biology of the investigation, rather than assay design.

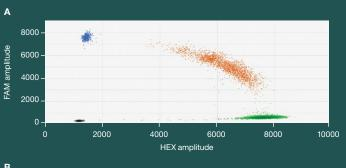
^{*} Each assay for the human, mouse, and rat genomes was experimentally tested for optimal efficiency, specificity, sensitivity, and linear dynamic range.

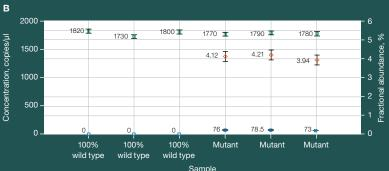
ABSOLUTE QUANTIFICATION USING EvaGreen



Droplet Digital PCR enables precise and reproducible detection calls across a range of sample concentrations. A, concentration plot shows data for both merged wells (solid blue markers) and individual replicate wells (open blue markers) across a dilution series. Interleukin 1 (*IL-1*) samples ranging from 0 to 50,000 copies were detected with an EvaGreen Assay and plotted on a log scale. There is excellent linearity and reproducibility between replicates. B, 1-D fluorescence amplitude plot shows a sample dilution curve detected with an EvaGreen *IL-1* Assay in the FAM/EvaGreen channel of the QX200 Droplet Reader. The data demonstrate good separation between positive () and negative () data points. All error bars are generated by QuantaSoft Software and represent a 95% confidence interval. NTC, no template control.

DETECTION OF EGFR L858R MUTATION





PrimePCR ddPCR Assays are fully validated for the detection of *EGFR L858* mutation. **A**, 2-D fluorescence amplitude plot shows triplicate wells of a mixed mutant and wild-type sample. The black cluster on the plot represents the negative droplets, the green cluster represents the droplets that are positive for wild-type DNA only, the blue cluster represents the droplets that are positive for mutant DNA only, and the orange cluster represents the droplets that are positive for both mutant and wild-type DNA. B, fractional abundance plot shows the percentage frequency (orange markers) of the mutant DNA in a wild-type DNA background. The blue markers indicate the concentration of mutant DNA (copies/µI) and the green markers indicate the concentration of wild-type DNA (copies/µl) in each of three replicate samples. All error bars are generated by QuantaSoft Software and represent a 95% confidence interval.



More Droplet Digital PCR Resources

Specifications			Catalog #	Description		
QX200 Droplet Generator Starting sample size 20 μl		1863021	200 x 20 μl reactions, 2x RT-ddPCR mix, includes			
Capacity Droplets per sa Dimensions (W	ample	1–8 samples/cartridge 20,000 28 x 36 x 13 cm (11 x 14 x 5 in.)	1863022	1 manganese acetate tube One-Step RT-ddPCR Kit for Probes, 5 ml (5 x 1 ml), 500 x 20 µl reactions, 2x RT-ddPCR mix, includes		
QX200 Drople Precision Linear dynamic Capacity Droplets per 96 Sample illumin	c range 6-well plate	±10% 5 orders of magnitude 1–96 samples Approximately 1,500,000 Light-emitting diodes	1864033 1864034 1864035	2 manganese acetate tubes QX200 ddPCR EvaGreen Supermix, 2 ml (2 x 1 ml), 200 x 20 µl reactions, 2x supermix QX200 ddPCR EvaGreen Supermix, 5 ml (5 x 1 ml), 500 x 20 µl reactions, 2x supermix QX200 ddPCR EvaGreen Supermix, 25 ml (5 x 5 ml),		
Sample detect Detection char Dimensions (W	ion nnels	Eight-enhithing diodes Multipixel photon counter FAM (EvaGreen), HEX (VIC) 66 x 52 x 29 cm (26 x 20 x 11 in.)	1864036	2,500 x 20 µl reactions, 2x supermix QX200 ddPCR EvaGreen Supermix, 50 ml (10 x 5 ml), 5,000 x 20 µl reactions, 2x supermix		
Ordering Information			1863052 1864052 1863040	ddPCR Buffer Control Kit, 9 ml (2 x 4.5 ml), 2x buffer QX200 Buffer Control for EvaGreen, 9 ml (2 x 4.5 ml), 2x buffer ddPCR Library Quantification Kit for Illumina TruSeq,		
Catalog # Description QX200 Droplet Digital PCR System 1864001 QX200 Droplet Digital PCR System, includes droplet generator, droplet reader, laptop computer, software, associated component		1863041	200 x 20 µl reactions, includes ddPCR Supermix for Probes (No dUTP) (2 x 1 ml vials), ddPCR Library Quantification Assay (1 x 200 µl vial), for quantification of Illumina TruSeq libraries using the QX200 or QX100 System ddPCR Library Quantification Kit for Ion Torrent,			
1864002	consumables QX200 Droplet Generator, includes droplet generator, 1 pkg of 24 DG8 Cartridges, 1 pkg of 24 DG8 Gaskets, 2 cartridge holders, 1 power cord QX200 Droplet Reader, includes droplet reader, ddPCR manual, 2 plate holders, USB cable, power cord Droplet Generator Cartridges and Gaskets, includes 5 pkg of 24 DG8 Cartridges, 5 pkg of 24 DG8 Gaskets DG8 Cartridges for QX200/QX100™ Droplet Generator, 1 pkg of 24 cartridges DG8 Gaskets for QX200/QX100 Droplet Generator, 1 pkg of 24 gaskets			200 x 20 µl reactions, includes ddPCR Supermix for Probes (No dUTP) (2 x 1 ml vials), ddPCR Library Quantification Assay (1 x 200 µl vial), for quantification of Ion Torrent libraries using the QX200 or QX100 System		
1864003			Thermal Cycler and Plate Sealer			
1864007			1851197	197 C1000 Touch™ Thermal Cycler with 96-Deep Well Reaction Module, includes C1000 Touch Thermal Cycler Chassis, 96-deep well reaction module, USB flash drive		
1864008			1814000	PX1 [™] PCR Plate Sealer, includes heat sealing instrument, plate support block that holds 96-well and 384-well plates, sealing frame, power cord		
1863009						
1863051 1863005	DG8 Cartridge Holder Droplet Generation Oil for Probes, 10 x 7 ml		EvaGreen is a trademark of Biotium, Inc. Bio-Rad Laboratories, Inc. is licensed by Biotium, Inc. to sell reagents containing EvaGreen Dye for use in real-time PCR,			

by Biotium, Inc. to sell reagents containing EvaGreen Dye for use in real-time PCR, for research purposes only. FAM, HEX, Ion Torrent, and VIC are trademarks of Life QX200 Droplet Generation Oil for EvaGreen, 2 x 7 ml Technologies Corporation. Illumina, MiSeg, and TruSeg are trademarks of Illumina, Inc. QX200 Droplet Generation Oil for EvaGreen, 10 x 7 ml Illumina is not affiliated with Bio-Rad Laboratories, Inc.

> The QX100 or QX200 Droplet Digital PCR System and/or its 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. Purchase of the product includes a limited, non-transferable right under such intellectual property for use of the product for internal research purposes only. No rights are granted for diagnostic uses. No rights are granted for use of the product for commercial applications of any kind, including but not limited to manufacturing, quality control, or commercial services, such as contract services or fee for services. 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.

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.

The QX200 Droplet Digital PCR System and Automated Droplet Generator are for research use only and are not intended for use in diagnostic procedures.





Bio-Rad Laboratories, Inc.

ddPCR Droplet Reader Oil, 2 x 1 L

2,500 x 20 µl reactions, 2x supermix

5,000 x 20 µl reactions, 2x supermix

200 x 20 µl reactions, 2x supermix

500 x 20 µl reactions, 2x supermix

2,500 x 20 µl reactions, 2x supermix

ddPCR Supermix for Probes, 25 ml (5 x 5 ml),

ddPCR Supermix for Probes, 50 ml (10 x 5 ml),

ddPCR Supermix for Probes (No dUTP), 2 ml (2 x 1 ml),

ddPCR Supermix for Probes (No dUTP), 5 ml (5 x 1 ml),

ddPCR Supermix for Probes (No dUTP), 25 ml (5 x 5 ml),

2x supermix

2x supermix

ddPCR Supermix for Probes, 2 ml (2 x 1 ml), 200 x 20 µl reactions,

ddPCR Supermix for Probes, 5 ml (5 x 1 ml), 500 x 20 µl reactions,

Life Science Group

1864005

1864006

1863004

1863026

1863010

1863027

1863028

1863023

1863024

1863025

ddPCR Reagents

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 Belgium 03 710 53 00 Brazil 55 11 3065 7550 Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 49 89 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300 Israel 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 64 9 415 2280 Norway 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 08 555 12700 Switzerland 026 674 55 05 Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000

Bulletin 6311 Rev C US/EG 15-0846 0615 Sig 1214