

# Counting DNA Molecule by Molecule

Adam S. Corner, PhD  
Bio-Rad Laboratories, Inc., Digital Biology Center, 5731 W Las Positas Blvd, Pleasanton, CA 94588



## Droplet Digital PCR

Bulletin 7374

### Abstract

Digital PCR is an established nucleic acid (NA) quantification technology, which enhances the broad use of PCR across molecular biology laboratories globally through improved precision and accuracy. Organizations that impact the quality of nucleic acid measurement and testing (metrology institutes, standards manufacturers, and quality assurance [QA]/proficiency testing organizations), adopted digital PCR early on and have since been pioneering its use for improved accuracy and precision in NA measurement. Ultimately, this will positively impact the quality of NA measurements generally and, more specifically, clinical NA testing. Molecular counting of nucleic acid molecules is the next measurement paradigm.

### Digital PCR — A Technology Step Change

Today digital PCR is a well-established technology, which enhances the broad use of PCR across molecular biology laboratories globally. From early attempts at limiting dilution PCR (Saiki et al. 1988) to using 384-well plates (Vogelstein and Kinzler 1999), through to microfluidic chambers on the Fluidigm Systems to current emulsion (Bio-Rad Laboratories, Inc., RainDance Technologies, Inc., Stilla Technologies) and chip-/plate-based approaches (Thermo Fisher Scientific Inc., JN Medsys, QIAGEN), digital PCR has become a variant of PCR adopted by those looking for greater precision and accuracy than that afforded by other amplification techniques, such as quantitative PCR (qPCR) and its variants. Among this range of platforms and formats, Bio-Rad has become the leader in the field with thousands of QX200 and QX ONE Droplet Digital PCR (ddPCR) Systems worldwide and more than 4,700 peer-reviewed publications covering applications spanning disease testing and monitoring, gene editing, cell therapies, and feed and food testing across the whole organismal spectrum from humans to viruses.

Often researchers progress from qPCR and similar amplification techniques to Droplet Digital PCR because of its technical benefits. For example, the use of Droplet Digital PCR in liquid biopsy analysis of circulating tumor DNA is realistically only possible at the current levels of sensitivity due to the partitioning of thousands of individual DNA molecules into thousands of individual droplets. Rare variant-containing molecules can then be detected in isolation within individual droplets. Rare variant detection that was previously restricted to >5% with qPCR and next-generation sequencing (NGS) is now readily achievable at <0.1% in a single well with highly sensitive Droplet Digital PCR.

Sensitivities of <0.02% (Lamy et al. 2015, Link-Lenczowska et al. 2018) for rare variants are possible within the population of 20,000 droplet partitions on the QX200 ddPCR System when utilizing single-well analyses. Levels of sensitivity can be further increased by merging multiple wells into a meta-well in an additive fashion (Hindson et al. 2011). Droplet Digital PCR thus has an expandable sensitivity range by adapting the number of wells utilized per sample, a definite benefit over qPCR or NGS.

**BIO-RAD**

Applications that utilize the precision of amplifying 20,000 droplet reactions in parallel have become core techniques for Droplet Digital PCR. Copy number variants that were typically restricted to determining haploid vs. diploid variants with qPCR can now precisely identify those ranging from 1 copy to >20 copies at the level of individual genome increments. Investigating copy number variants associated with cancer development and progression, for example, *HER2* monitoring in breast cancer, can be challenging and open to subjective interpretation. With Droplet Digital PCR, clinical interpretation of *HER2* amplification between <2 copies vs. >2 copies is readily achievable (Garcia-Murillas et al. 2013, Suryavanshi et al. 2019).

Similarly, ascertaining copy number variants beyond haploid vs. diploid variants in germ line scenarios is achievable, for example, distinguishing *SMN1* and *SMN2* gene copy variations in newborns diagnosed with spinal muscular atrophy (Vidal-Folch et al. 2018, Chien et al. 2017) and identifying gene copy variation as a contributing factor in schizophrenia development (Marshall et al. 2017).

In additional research, the effort to quantify copy number variations (CNVs) in polyploidy plant species (Juoanin et al. 2020) and determine transgene copy number in plants (Collier et al. 2017) benefits greatly from the precision achievable with Droplet Digital PCR.

Ultimately bringing together both the sensitivity and precision of Droplet Digital PCR, CNV detection in rare genomic fractions is possible with noninvasive prenatal testing and graft vs. host disease detection in patients with transplants.

Detecting small amounts of fetal DNA is technically challenging and yet achievable within the large background of maternal DNA (typically 4–5% fetal) using Droplet Digital PCR. But being able to detect small changes (2 vs. 3 chromosome copies for fetal aneuploidy) within the fetal subpopulation of maternal DNA requires combined levels of precision and sensitivity unachievable with other PCR techniques. This application does require >20,000 droplets. But with the meta-well merging capabilities of Droplet Digital PCR described above, detection of fetal aneuploidy is readily achievable (El Khattabi et al. 2016, Lee and Hwang 2015).

Similarly, several different methods have been described to detect donor-derived circulating tumor DNA (ddctDNA) in the blood of recipients with transplants using Droplet Digital PCR. Goh et al. (2017) utilized donor vs. recipient discriminatory insertion-deletion mutations, Zou et al. (2017) utilized human leukocyte antigen-specific probe sets to identify ddctDNA, and Oellerich et al. (2014) utilized a comprehensive panel of single nucleotide polymorphisms to identify ddctDNA in recipient plasma. These methods again utilize the ddPCR

technical benefits of increased precision and sensitivity over alternative PCR-based approaches to allow for the monitoring of small amounts of circulating DNA.

More recently, determination of viral copy number (VCN) of chimeric antigen receptor T (CAR-T) cells and T-cell receptor (TCR) cell lines by Droplet Digital PCR has been published on extensively (Fehse et al. 2020, Kao et al. 2018, Lu et al. 2020). Precision monitoring of adeno-associated virus and lentiviral integration into cells for downstream gene therapy use in human clinical treatments is a huge regulatory and clinical issue, as measuring VCN accurately is critical in producing the optimal dose for patients. From both technical and clinical perspectives there is growing acceptance that Droplet Digital PCR offers a fully viable route to determining VCN for clinical dosing of CAR-T cells and TCR cells within recommended safe ranges due to its absolute quantification capabilities and precision.

### **The Science of Measurement**

In parallel with the research groups using Droplet Digital PCR to further our understanding of biology across the globe and enhance our clinical treatments of the human condition, there is a research community that underpins every measurement that is made in every laboratory. Metrology is the science of measurement and forms the foundation of the vast majority of processes that occur throughout a normal day, for example, the speed and distance you travel to work, the volume of coffee you buy on the way to work, and the weight and electrical power of the elevator that takes you up to your office! The majority of the units of measurement are founded on Système International (SI) units governed by the global measurement organization, the Bureau International des Poids et Mesures. Implementation of these units of measurement is monitored, managed, and promoted by metrology laboratories within each country.

In molecular biology and clinical and research studies, metrology labs were early adopters and fully implemented digital PCR with a focus on improving precision and accuracy in measuring nucleic acids. The biological application goal across global metrology labs, for example, for liquid biopsy variants, viral detection, and genetically modified organism (GMO) detection, is almost secondary to the more global goal of defining good measurement practices and the development of good reference materials. Each National Institute of Metrology tends to have a different biological focus, or several, but a common metrological focus to improve measurement processes and practices in science.

The National Measurement Institute of Australia has focused on the applications of nucleic acid copy number count, copy number variation, and absolute nucleic acid quantification, key techniques for which Droplet Digital PCR is well-suited. From the very earliest of papers on the use of Droplet Digital PCR

to measure lambda DNA quantities (Pinheiro et al. 2012) and assessment of droplet volumes, in collaboration with the Joint Research Center of the European Commission, to improve absolute quantification with Droplet Digital PCR (Corbisier et al. 2015) to more recent work on the development of a reference material for quality control of Droplet Digital PCR for absolute nucleic acid quantification (Pinheiro et al. 2017), Australian metrology studies have confirmed it as a suitable technology for absolute quantification of nucleic acids.

The UK National Measurement Laboratory has published on a diverse range of applications, covering liquid biopsy *KRAS* variants (Whale et al. 2016), influenza and human immunodeficiency virus (HIV) detection (Whale et al. 2016), and bacterial detection (Devonshire et al. 2015), often with a clinical focus on the application of measurement. The work at the National Measurement Laboratory also highlighted the robust repeatability and reproducibility of Droplet Digital PCR as well as its ability to perform absolute quantification without standards or reference materials. Raising awareness of the inherent challenges of using standards for qPCR was particularly well illustrated by the work on the HIV calibration standard 8E5 (Busby et al. 2017).

The National Institute of Biology in Slovenia has published several articles about the detection and quantification of genetically modified organisms (Dobnik et al. 2015, Kosir et al. 2019) and viral detection (Dobnik et al. 2019, Mehle et al. 2018, Pavsic et al. 2017). This work highlighted that Droplet Digital PCR, a relatively new technology, can be heavily multiplexed with no loss of precision or accuracy compared to qPCR. It also demonstrated (with other metrology institutes) that Droplet Digital PCR has high reproducibility and repeatability, without the use of a standard curve, and, as such, is potentially suitable as a metrological reference method. These two factors (multiplexing and lack of standards) are significant for future laboratories' adoption of Droplet Digital PCR in place of established qPCR. This latter point is well illustrated by the relatively early adoption of Droplet Digital PCR by groups aiming to determine VCN in the gene editing field.

In support of the use of Droplet Digital PCR for reference material validation, the National Institute of Standards and Technology (NIST) in the U.S. has utilized digital PCR to carry out extensive research on absolute quantification of both double- and single-stranded DNA (Dewer et al. 2018, Kline and Dewer 2017, Kline and Dewer 2019). Interestingly, although much of the work around Droplet Digital PCR has been focused on defining the technology as being suitable for traceable nucleic acid quantification, NIST also

has been working with digital PCR in various forms for the development of reference standards over the last 8 years. Early microchamber digital PCR systems were utilized to develop viral reference materials for cytomegalovirus (Haynes 2013). More recently, ddPCR technologies were adopted to develop further reference materials for the BK virus (Cleveland et al. 2018) as well as in oncology for *HER2* gene amplification detection and *EGFR* and *MET* gene copy number analysis (He et al. 2016, He et al. 2019), and in the development of an international reference material for quantification of human genomic DNA (Romsos et al. 2018).

Similarly to NIST in the U.S., the Joint Research Center of the European Commission has been involved in metrological determination of reference materials and nucleic acid practices across a range of applications, including validation of GMO control plasmid for qPCR testing (Grohmann et al. 2017), copy number determination (Corbisier et al. 2015, Deprez et al. 2016), and liquid biopsy ctDNA analysis (Whale et al. 2018). But very early on in their adoption of digital PCR, these researchers highlighted the potential of the technique for analytical assessment of nucleic acids.

In the last 8 years, early publications by metrology labs have focused on assessing digital PCR for accurate and precise measurement of nucleic acids by comparing it to qPCR. Quantitative PCR has the potential for high precision but has limitations around accuracy and a large potential for bias (uncertainty) when using poor-quality standards for a standard curve. Droplet Digital PCR has become established as a technique with the precision to match qPCR and absolute accuracy better than that of qPCR. It provides absolute quantification without the need for standards.

Metrology experiments are often structured around high levels of intralab and interlab replicates to illustrate that Droplet Digital PCR is robust for comparison within and between labs, again without the need for common or international standards. Test samples are typically set up gravimetrically with the aim of reducing experimental uncertainty, thus ensuring accuracy, which leads to a better understanding of the underlying repeatability and reproducibility of the ddPCR technique. Finally, comparison to orthogonal technologies for measuring nucleic acids has taken Droplet Digital PCR closer to being defined as an independent, traceable technique for nucleic acid quantification (Whale et al. 2018, Yoo et al. 2016). This idea was also proposed by Mehle et al. (2018) with Droplet Digital PCR compared to qPCR and Kline and Dewer (2017) in their comparison of chip-based digital PCR (cdPCR) to Droplet Digital PCR.

## Molecular Counting

By applying Droplet Digital PCR to a range of real-world application challenges, the National Measurement Laboratories in multiple countries have illustrated that digital PCR has the following properties:

- **Repeatability** — intralab studies illustrate the same answer, without the need for calibrators or standard curves
- **Reproducibility** — interlab studies illustrate the same answer, without standards or calibrators, or even the same assay design
- **Reduced uncertainty (absolute quantification)** — by comparison to orthogonal methods (for example, inductively coupled plasma–mass spectrometry and cdPCR), the molecular count answer is as close to truth as possible with minimal (and acceptable) uncertainty

And as an outcome of these technological capabilities Droplet Digital PCR has been accepted as a primary reference measurement procedure for SI traceable quantification of nucleic acid primary reference materials through molecular enumeration (ISO 20395:2019, ISO 17511:2020). By comparison, qPCR is recognized as a traceable method, but relies on suitable traceable reference materials.

Droplet Digital PCR can be used to simply count your DNA molecules (taking into consideration specificity, uncertainty, and completeness of count)!

### Manufacturers of Reference and Quality Control Standards

Besides the metrology labs aiming to define good metrological practices in molecular biology and rapidly adopting Droplet Digital PCR for that purpose in nucleic acid quantification, there are organizations that are utilizing the benefits of Droplet Digital PCR at a more applied level.

In the United Kingdom, the National Institute for Biological Standards and Controls (NIBSC) plays a major role in production and distribution of international reference materials, including the majority of World Health Organization (WHO) international standards and reference materials. In the context of nucleic acid detection and testing, NIBSC produces a wide range of viral and bacterial WHO controls that are defined based on the International Unit, an accepted arbitrary unit of definition. These controls are typically defined by a large interlab study using an extensive range of nucleic acid detection techniques to illustrate broad commutability of the control (NIBSC code 09/162; Fryer et al. 2016). NIBSC also produces genomic controls utilizing Droplet Digital PCR as

a key technology during validation in recognition of its inherent precision and absolute quantitative capability for nucleic acids (NIBSC code 18/130: WHO International Standard, Sanzone et al. 2019).

Manufacturers of quality control (QC) standards have been relatively early adopters of Droplet Digital PCR, particularly in the areas of translational research and oncology. One of the major technical benefits of Droplet Digital PCR is the ability to precisely measure the proportion of a rare DNA sequence in the presence of a common variant. This is perfectly exemplified when monitoring somatic variants in the blood of patients with cancer. ctDNA can be present in low-integer numbers in a background of thousands of wild-type DNA molecules. Droplet Digital PCR is well adapted for detecting these rare fragments.

However, to validate this process in translational and clinical laboratories, controls must be utilized under current testing paradigms to validate the sensitivity of the ddPCR platform. To that end, companies such as SeraCare Life Sciences, SensiD GmbH, and Horizon Discovery Ltd. have developed a range of oncology controls with mutant fractions applicable across a broad range (<0.1–10%) for this purpose and use Droplet Digital PCR as their QC technology. Exact Diagnostics, which produces control materials for the virology and bacteriology markets, utilizes Droplet Digital PCR for its precision and accuracy in the infectious disease field, where absolute accuracy and sensitivity are critical requirements.

### External Quality Assurance and Proficiency Testing

Moving closer toward the clinical testing world, there is a recognition among external quality assurance (EQA) and proficiency testing (PT) organizations that Droplet Digital PCR may offer benefits for nucleic acid detection and quantification. This may come from the indirect benefit of buying EQA controls from an established QC manufacturer that has adopted Droplet Digital PCR (for example, SeraCare *EGFR* Panel for Genomics Quality Assessment [GenQA]) and, thus, providing higher-quality reference materials for an EQA. Or it may come from the direct adoption of Droplet Digital PCR into the development path of an EQA either during the production of internal EQA controls (for example, Quality Control for Molecular Diagnostics [QCMD]) or in the quantification of participant samples as part of an NA extraction EQA (GenQA). All of these approaches improve accuracy and precision across the EQA process.

The purpose of EQA organizations is not to promote individual technologies, but to assess the use of a range of technologies used by clinical labs and support those labs in achieving suitable QA accreditation. But recognition of the benefits of digital PCR over other nucleic acid amplification test technologies may present itself through the cumulative results generated over future EQA schemes. Further recognition of this benefit by participating clinical labs will ultimately result in better and safer clinical testing.

### **The Next Steps for Absolute Nucleic Acid Quantification**

Absolute quantification by qPCR requires standards to generate a standard curve, of known quantity, which, ideally, is SI traceable. Variant allele frequencies determined by NGS require controls and calibrators that confirm mutant allele ratios and limits of detection and quantification. Since there is a large establishment of testing labs equipped with qPCR systems and NGS systems, accurate standards generated by Droplet Digital PCR do reduce overall measurement uncertainty in an SI traceable manner. While pragmatically accepting that many labs cannot discontinue old technology overnight for the next-generation technology, this approach does endorse Droplet Digital PCR as the core NA quantifying technology that subsequently supports the qPCR and NGS instrument/testing base.

Over time, many core labs may choose to adopt Droplet Digital PCR directly rather than obtain SI traceable standards for use as controls for second-generation qPCR technologies. There is already growing literature illustrating NGS labs utilizing Droplet Digital PCR as not only an orthogonal validation method but also indirectly as a measurement technique that is SI traceable.

### **Conclusions**

Although a relatively unknown sector within molecular biology and clinical testing, it is vitally important that metrologists (who are molecular biologists) both validate new technologies for NA quantification and promote the benefits of those technologies if proven. It is relatively easy to understand that quantifying DNA (or RNA) using Droplet Digital PCR may improve downstream manipulation of samples in subsequent molecular biology experiments, for example, input sample

amounts for NGS library preparation. However, the more important purpose lies in understanding how being able to quantify NA more accurately, as defined by National Measurement Laboratories, has a direct impact on standards manufacturers and EQA and PT organizations, and, ultimately, on clinical labs that are diagnosing human disease.

As Droplet Digital PCR becomes established as an independent reference measurement technology, its value as an orthogonal (SI traceable) validation technology for NGS also becomes recognized. This, in turn, impacts the quality of clinical testing by NGS panels. Beyond high value clinical testing (NGS validation and quantification of nucleic acids), where the benefits of adoption support the added costs of the technology over qPCR, adoption by clinical and research labs on a broader scale must balance the costs of adopting new technology against the benefits of improved precision and absolute quantification. Ultimately, there are opportunities for digital PCR to become a routine NA quantification method, in a format that is smaller, less expensive, and with a faster turnaround time than today.

There is recognition within the scientific community as a whole that Droplet Digital PCR offers improved accuracy and precision over qPCR, and this leads to the ability to determine low-level variant allele frequencies and precise CNV detection. But despite broad recognition of Droplet Digital PCR for independent absolute quantification, adoption is still not widespread, even though the majority of publications illustrating Droplet Digital PCR cite “absolute quantification without standards” as a major benefit over other NA amplification technologies in their introductions. This reticence to count molecules absolutely is illustrated clearly by the majority of translational research publications in oncology investigating somatic mutations and declaring the mutant copy number as a percentage or ratio to wild-type genomes rather than mutant copies per unit of blood or plasma. This may change if Droplet Digital PCR is accepted as a primary reference measurement procedure for SI traceable quantification of primary reference materials through molecular enumeration. This would open the door for Droplet Digital PCR to become a primary reference measurement procedure in the routine laboratory by molecular counting.

## References

- Busby E et al. (2017). Instability of 8E5 calibration standard revealed by digital PCR risks inaccurate quantification of HIV DNA in clinical samples by qPCR. *Sci Rep* 7, 1,209.
- Chien Y-H et al. (2017). Presymptomatic diagnosis of spinal muscular atrophy through newborn screening. *J Pediatr* 190, 124–129.
- Cleveland MH et al. (2018). Certification of Standard Reference Material 2365 BK Virus DNA Quantitative Standard. National Institute of Standards and Technology. U.S. Department of Commerce. NIST Standard Reference Materials Publication 260-191.
- Collier R et al. (2017). Accurate measurement of transgene copy number in crop plants using droplet digital PCR. *Plant J* 90, 1,014–1,025.
- Corbisier P et al. (2015). DNA copy number concentration measured by digital and droplet digital quantitative PCR using certified reference materials. *Anal Bioanal Chem* 407, 1,831–1,840.
- Deprez L et al. (2016). Validation of a digital PCR method for quantification of DNA copy number concentrations by using a certified reference material. *Biomol Detect Quantif* 30, 29–39.
- Devonshire AS et al. (2015). Highly reproducible absolute quantification of *Mycobacterium tuberculosis* complex by digital PCR. *Anal Chem* 87, 3,706–3,713.
- Dobnik D et al. (2015). Multiplex quantification of 12 European Union authorized genetically modified maize lines with droplet digital polymerase chain reaction. *Anal Chem* 87, 8,218–8,226.
- Dobnik D et al. (2019). Accurate quantification and characterization of adeno-associated viral vectors. *Front Microbiol* 10, 1,570.
- Duewer DL et al. (2018). Evaluating droplet digital PCR for the quantification of human genomic DNA: Converting copies per nanoliter to nanograms nuclear DNA per microliter. *Anal Bioanal Chem* 410, 2,879–2,887.
- El Khattabi LA et al. (2016). Could digital PCR be an alternative as a non-invasive prenatal test for trisomy 21: A proof of concept study. *PLoS One* 11, e0155009.
- Fehse B et al. (2020). Digital PCR assays for precise quantification of CD19-CAR-T cells after treatment with axicabtagene ciloleucel. *Mol Ther Methods Clin Dev* 16, 172–178.
- Fryer JF et al. (2016). A collaborative study to establish the 1st WHO International Standard for human cytomegalovirus for nucleic acid amplification technology. *Biologicals* 44, 242–251.
- Garcia-Murillas I et al. (2013). Determination of HER2 amplification status on tumour DNA by digital PCR. *PLoS One* 8, e83409.
- Goh SK et al. (2017). Probe-free digital PCR quantitative methodology to measure donor-specific cell-free DNA after solid-organ transplantation. *Clin Chem* 63, 742–750.
- Grohmann L et al. (2017). Screening for six genetically modified soybean lines by an event-specific multiplex PCR method: Collaborative trial validation of a novel approach for GMO detection. *J Consum Prot Food Saf* 12, 23–36.
- Haynes RJ et al. (2013). Standard reference material 2366 for measurement of human cytomegalovirus DNA. *J Mol Diagn* 15, 177–185.
- He H-J et al. (2016). Development of NIST standard reference material 2373: Genomic DNA standards for HER2 measurements. *Biomol Detect Quantif* 8, 1–8.
- He H-J et al. (2019). Development and interlaboratory evaluation of a NIST reference material RM 8366 for EGFR and MET gene copy number measurements. *Clin Chem Lab Med* 57, 1,142–1,152.
- Hindson BJ et al. (2011). High-throughput droplet digital PCR system for absolute quantitation of DNA copy number. *Anal Chem* 83, 8,604–8,610.
- ISO 17511:2020. In vitro diagnostic medical devices — Requirements for establishing metrological traceability of values assigned to calibrators, trueness control materials and human samples.
- ISO 20395:2019. Biotechnology — Requirements for evaluating the performance of quantification methods for nucleic acid target sequences — qPCR and dPCR.
- Juoanin A et al. (2020). Optimisation of droplet digital PCR for determining copy number variation of  $\alpha$ -gliadin genes in mutant and gene-edited polyploid bread wheat. *J Cereal Sci* 92, 102903.
- Kao RL et al. (2018). A cetuximab-mediated suicide system in chimeric antigen receptor-modified hematopoietic stem cells for cancer therapy. *Hum Gene Ther* 30, 413–428.
- Kline MC and Duewer DL (2017). Evaluating droplet digital polymerase chain reaction for the quantification of human genomic DNA: Lifting the traceability fog. *Anal Chem* 89, 4,648–4,654.
- Kline MC and Duewer DL (2019). NIST Special Publication 1200-27: Evaluation of Methods for Assessing the Proportion of Single Stranded Nuclear DNA in Human Blood Extracts. National Institute of Standards and Technology. U.S. Department of Commerce.
- Kosir AB et al. (2019). Digital PCR as an effective tool for GMO quantification in complex matrices. *Food Chem* 294, 73–78.
- Lamy P-J et al. (2015). Next-generation genotyping by digital PCR to detect and quantify the BRAF V600E mutation in melanoma biopsies. *J Mol Diagn* 17, 366–373.
- Lee SY and Hwang SY (2015). Application of digital polymerase chain reaction technology for noninvasive prenatal test. *J Genet Med* 12, 72–78.
- Link-Lenczowska D et al. (2018). A comparison of qPCR and ddPCR used for quantification of the JAK2 V617F allele burden in Ph negative MPNs. *Ann Hematol* 97, 2,299–2,308.
- Lu A et al. (2020). Application of droplet digital PCR for the detection of vector copy number in clinical CAR/TCR T cell products. *J Transl Med* 18, 191.
- Marshall CR et al. (2017). Contribution of copy number variants to schizophrenia from a genome-wide study of 41,321 subjects. *Nat Genet* 49, 27–35.
- Mehle N et al. (2018). Validated reverse transcription droplet digital PCR serves as a higher order method for absolute quantification of Potato virus Y strains. *Anal Bioanal Chem* 410, 3,815–3,825.
- Oellerich M et al. (2014). Use of graft-derived cell-free DNA as an organ integrity biomarker to reexamine effective tacrolimus trough concentrations after liver transplantation. *Ther Drug Monit* 36, 136–140.
- Pavsic J et al. (2017). Inter-laboratory assessment of different digital PCR platforms for quantification of human cytomegalovirus DNA. *Anal Bioanal Chem* 409, 2,601–2,614.
- Pinheiro LB et al. (2012). Evaluation of a droplet digital polymerase chain reaction format for DNA copy number quantification. *Anal Chem* 84, 1,003–1,011.
- Pinheiro LB et al. (2017). Interlaboratory reproducibility of droplet digital polymerase chain reaction using a new DNA reference material format. *Anal Chem* 89, 11,243–11,251.
- Romsos EL et al. (2018). Certification of Standard Reference Material 2372a Human DNA Quantification Standard. National Institute of Standards and Technology. U.S. Department of Commerce. NIST Standard Reference Materials Publication 260-189.
- Saiki RK et al. (1988). Primer directed enzymatic amplification of DNA with a thermostable DNA polymerase. *Science* 239, 487–491.
- Sanzone AP et al. (2019). Collaborative Study to Evaluate the Proposed WHO 1st International Standards for Cancer Genomes. World Health Organization. WHO/BS/2019.2368.
- Suryavanshi M et al. (2019). Droplet digital polymerase chain reaction offers an improvisation over conventional immunohistochemistry and fluorescent *in situ* hybridization for ascertaining Her2 status of breast cancer. *South Asian J Cancer* 8, 203–210.

Vidal-Folch N et al. (2018). Multiplex droplet digital PCR method applicable to newborn screening, carrier status, and assessment of spinal muscular atrophy. *Clin Chem* 64, 1,753–1,761.

Vogelstein B and Kinzler KW (1999). Digital PCR. *Proc Natl Acad Sci USA* 96, 9,236–9,241.

Whale AS et al. (2016). Detection of rare drug resistance mutations by digital PCR in a human influenza A virus model system and clinical samples. *J Clin Microbiol* 54, 392–400.

Whale AS et al. (2016). International interlaboratory digital PCR study demonstrating high reproducibility for the measurement of a rare sequence variant. *Anal Chem* 89, 1,724–1,733.

Whale AS et al. (2018). Assessment of digital PCR as a primary reference measurement procedure to support advances in precision medicine. *Clin Chem* 64, 1,296–1,307.

WHO International Standard: 1st International Standard 2019 for MOLT-4 Cancer Genome. NIBSC code 18/130. Version 1.0. March 16, 2020. National Institute for Biological Standards and Control. World Health Organization.

Yoo H-B et al. (2016). International comparison of enumeration-based quantification of DNA copy-concentration using flow cytometric counting and digital polymerase chain reaction. *Anal Chem* 88, 12,169–12,176.

Zou J et al. (2017). Rapid detection of donor cell free DNA in lung transplant recipients with rejections using donor-recipient HLA mismatch. *Hum Immunol* 78, 342–349.

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

BIO-RAD, DDPCR, DROPLET DIGITAL PCR, and QX ONE are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

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 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](https://www.bio-rad.com/en-us/trademarks) for details.

All trademarks used herein are the property of their respective owner.



**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

