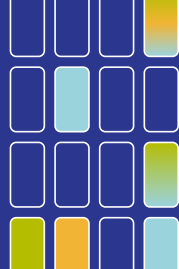


Comparing real-time and digital PCR technologies

Both digital PCR (dPCR) and real-time or quantitative PCR (qPCR) can be used to quantify nucleic acids in a sample. This is performed by amplifying a target nucleic acid molecule with a DNA polymerase enzyme.

dPCR

Distribute reaction, amplify, and count at endpoint

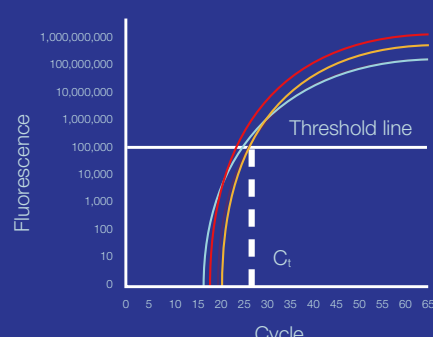


Absolute measurement—counts target of interest via single-molecule amplification across a large number of PCR replicates. Run at limiting dilution to ensure at least one reaction does not contain target DNA.

vs.

qPCR

Measure bulk reaction fluorescence at each cycle until plateau phase



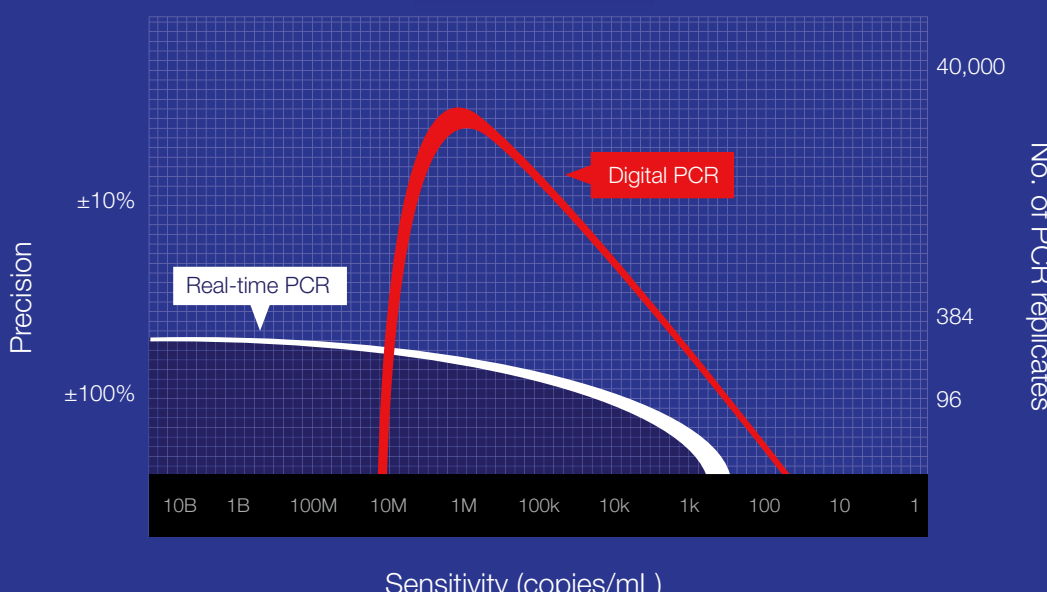
Measures PCR amplification against a reference as it occurs. Data are collected during the exponential (log) phase of PCR.



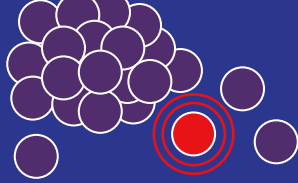
TECHNOLOGY



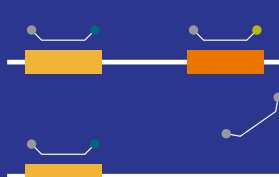
QUANTITATIVE



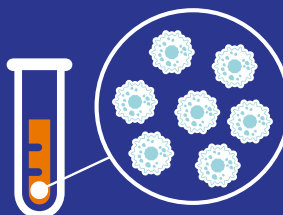
APPLICATIONS



Rare-target detection



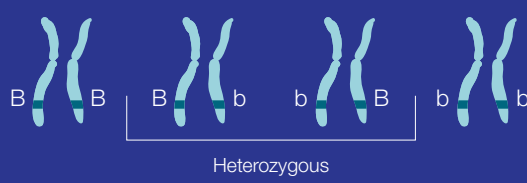
Single molecule characterization and quantification



Absolute quantification of viral load



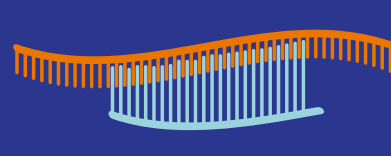
Somatic copy number variation or low fold changes



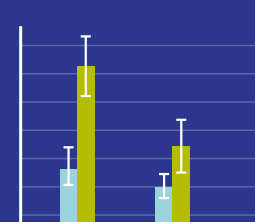
Heterozygous
SNP genotyping



Relative gene expression analysis



MicroRNA analysis



Standard copy number variation



WORKFLOW

Mix PCR reagents: nucleic acid, master mix, primers, and probes



Load wells of plate



Place on instrument for dPCR

[Watch the video](#)

Mix PCR reagents: nucleic acid, master mix, primers, and probes



Load individual strip tubes or plate



Place on instrument for qPCR



ADVANTAGES

Quantitative data output—no reliance on references or standards for conversion of data points

Capable of analyzing rare targets against wild-type or non-target background

Unlike traditional qPCR, digital PCR provides a linear response to the number of copies present to allow for small fold-change differences to be detected

Single molecule resolution interrogation enables identification and quantification of molecules containing multiple targets (e.g., phased targets or engineered plasmids)

Improved tolerance to some PCR inhibitors

Broadly accepted, well-established protocols and assays

Increased dynamic range of detection

Detection is capable down to a 2-fold change

Higher sample throughput with lower cost

Collects data in the exponential phase of PCR, providing a permanent record of amplicon amplification