



qPCR and dPCR for biopharma

## QualTrak real-time PCR and digital PCR ecosystem for streamlined biologics development

Biologics—including vaccines, cell and gene therapies, monoclonal antibodies (mAbs), and their biosimilars—have great potential for the treatment and prophylaxis of a broad range of diseases. As of 2018, eight of the top ten best-selling drugs in the US were mAbs [1]. mAb biosimilars are expected to follow a comparable path, with the biosimilars market set to double by 2025 [2]. Cell and gene therapies are also predicted to increase, with a projected compound annual growth rate (CAGR) of almost 25% in the near future [3]. This growth offers tremendous opportunities, especially for manufacturers who can develop new and effective therapies quickly.

Rapid and efficient development of new and effective biologics is no easy feat. Researching, developing, and manufacturing new biologic therapies is an arduous and expensive process fraught with regulatory and supply chain challenges. It takes an average of 10 years and \$2.6 billion to develop a new treatment, and only 12% of candidates make it to market [4], while others end up in “the valley of death”—the transition from laboratory to trial where candidates often fail.

With such drastic implications for time and resources, it is imperative that biologic manufacturers choose the right commercial collaborators from the start—agile partners who can improve predictivity, speed, and consistency to bridge basic and clinical research.

## Work with a partner who offers reliable, high-quality PCR products for every step of the development process

### The role of Thermo Fisher Scientific in PCR-based detection of coronavirus

“PCR” became a household term because of its common use in SARS-CoV-2 detection and surveillance; however, its applications in the coronavirus crisis do not end there. In particular, qPCR continues to play an important role in the development of SARS-CoV-2 treatments and vaccines. qPCR assays, reagents, and instruments developed by Thermo Fisher were not only cited in detection, but also in the development research for several SARS-CoV-2 vaccines, contributing on many levels to pandemic mitigation efforts.

### Leverage high-throughput, high-quality products throughout the entire workflow

#### Thanks for reading!

Access the unabridged PDF and more at the QualTrak resources page.

Keep reading



Table 1. Suggested applications for qPCR and dPCR in biopharma discovery, development, and manufacturing.

Phase	Goal	mAbs applications	Vaccine applications	Cell and gene therapy applications
Discovery	<ul style="list-style-type: none"> <li>Target identification</li> <li>Target validation</li> <li>Target characterization</li> </ul>	<ul style="list-style-type: none"> <li>Target identification</li> <li>Target validation</li> <li>Target characterization</li> </ul>	<ul style="list-style-type: none"> <li>Target identification</li> <li>Target validation</li> <li>Target characterization</li> </ul>	<ul style="list-style-type: none"> <li>Target identification</li> <li>Target validation</li> <li>Target characterization</li> </ul>
Process development	<ul style="list-style-type: none"> <li>Process optimization</li> <li>Scale-up</li> <li>Quality control</li> </ul>	<ul style="list-style-type: none"> <li>Process optimization</li> <li>Scale-up</li> <li>Quality control</li> </ul>	<ul style="list-style-type: none"> <li>Process optimization</li> <li>Scale-up</li> <li>Quality control</li> </ul>	<ul style="list-style-type: none"> <li>Process optimization</li> <li>Scale-up</li> <li>Quality control</li> </ul>
Preclinical and clinical research	<ul style="list-style-type: none"> <li>Target validation</li> <li>Target characterization</li> <li>Target identification</li> </ul>	<ul style="list-style-type: none"> <li>Target validation</li> <li>Target characterization</li> <li>Target identification</li> </ul>	<ul style="list-style-type: none"> <li>Target validation</li> <li>Target characterization</li> <li>Target identification</li> </ul>	<ul style="list-style-type: none"> <li>Target validation</li> <li>Target characterization</li> <li>Target identification</li> </ul>
Manufacturing	<ul style="list-style-type: none"> <li>Quality control</li> <li>Scale-up</li> <li>Process optimization</li> </ul>	<ul style="list-style-type: none"> <li>Quality control</li> <li>Scale-up</li> <li>Process optimization</li> </ul>	<ul style="list-style-type: none"> <li>Quality control</li> <li>Scale-up</li> <li>Process optimization</li> </ul>	<ul style="list-style-type: none"> <li>Quality control</li> <li>Scale-up</li> <li>Process optimization</li> </ul>

**Thanks for reading!**  
 Access the unabridged PDF and more at the QualTrack resources page.

[Keep reading](#)

# qPCR and dPCR offer more consistent results due to innovative instrumentation, optimized master mixes, and predesigned and custom assays

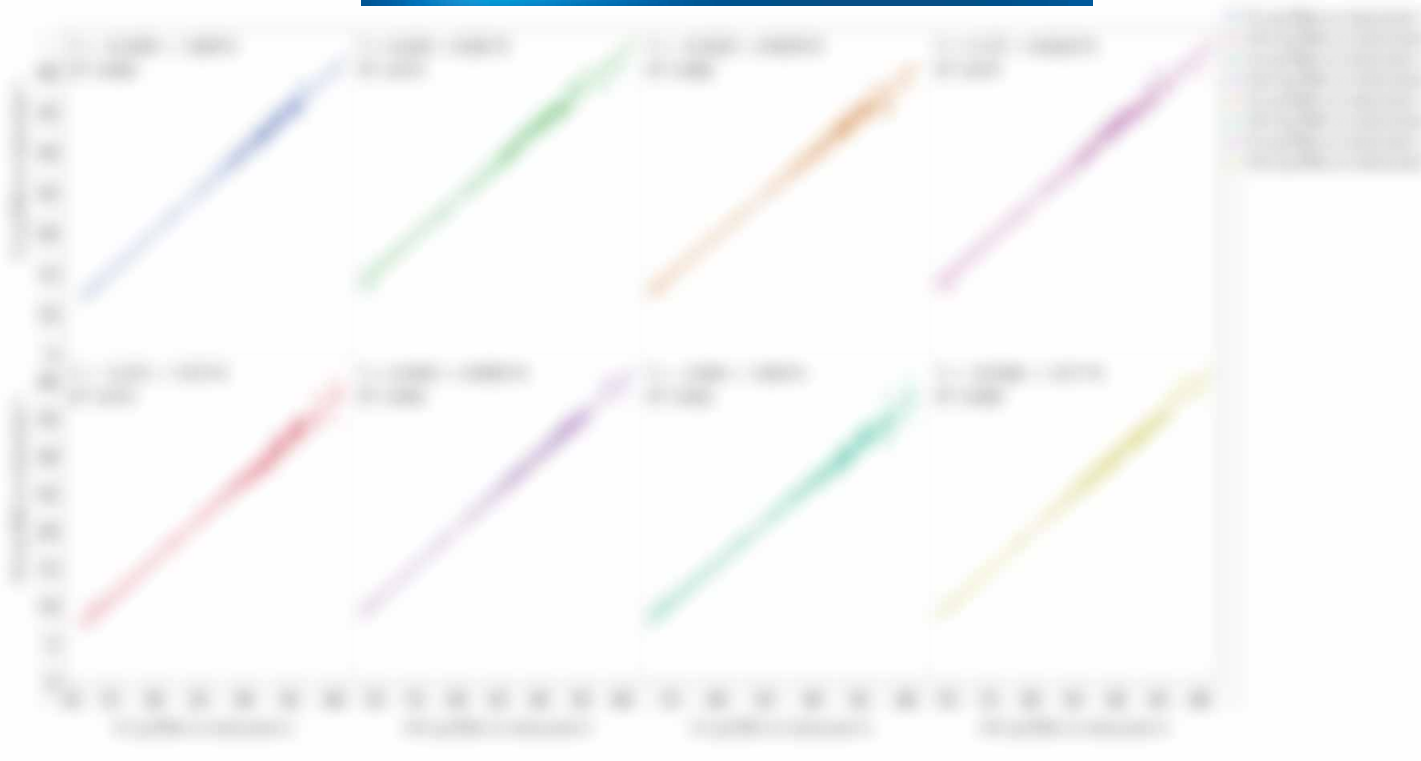
## Instrumentation

The instrumentation used in qPCR and dPCR is highly sensitive and accurate, allowing for the detection of low levels of target DNA. This is achieved through the use of highly sensitive detection systems, such as fluorescence and electrochemiluminescence. The instrumentation is also highly precise, allowing for the detection of small differences in target DNA levels. This is achieved through the use of highly precise amplification and detection systems, such as digital PCR and microfluidics. The instrumentation is also highly flexible, allowing for the detection of a wide range of target DNA sequences. This is achieved through the use of highly flexible amplification and detection systems, such as multiplexed qPCR and digital PCR.

The instrumentation used in qPCR and dPCR is highly sensitive and accurate, allowing for the detection of low levels of target DNA. This is achieved through the use of highly sensitive detection systems, such as fluorescence and electrochemiluminescence. The instrumentation is also highly precise, allowing for the detection of small differences in target DNA levels. This is achieved through the use of highly precise amplification and detection systems, such as digital PCR and microfluidics. The instrumentation is also highly flexible, allowing for the detection of a wide range of target DNA sequences. This is achieved through the use of highly flexible amplification and detection systems, such as multiplexed qPCR and digital PCR.

**Thanks for reading!**  
Access the unabridged PDF and more at the QualTrack resources page.

[Keep reading](#)



**Figure 1. Reproducibility of gene expression measurements using Applied Biosystems™ MicroAmp™ Fast Optical 96-Well Reaction Plates on different instruments.** Pairwise comparisons of cDNA data from 10 ng or 100 ng universal human reference (UHR) RNA on 3 different instruments show high correlations from instrument to instrument.



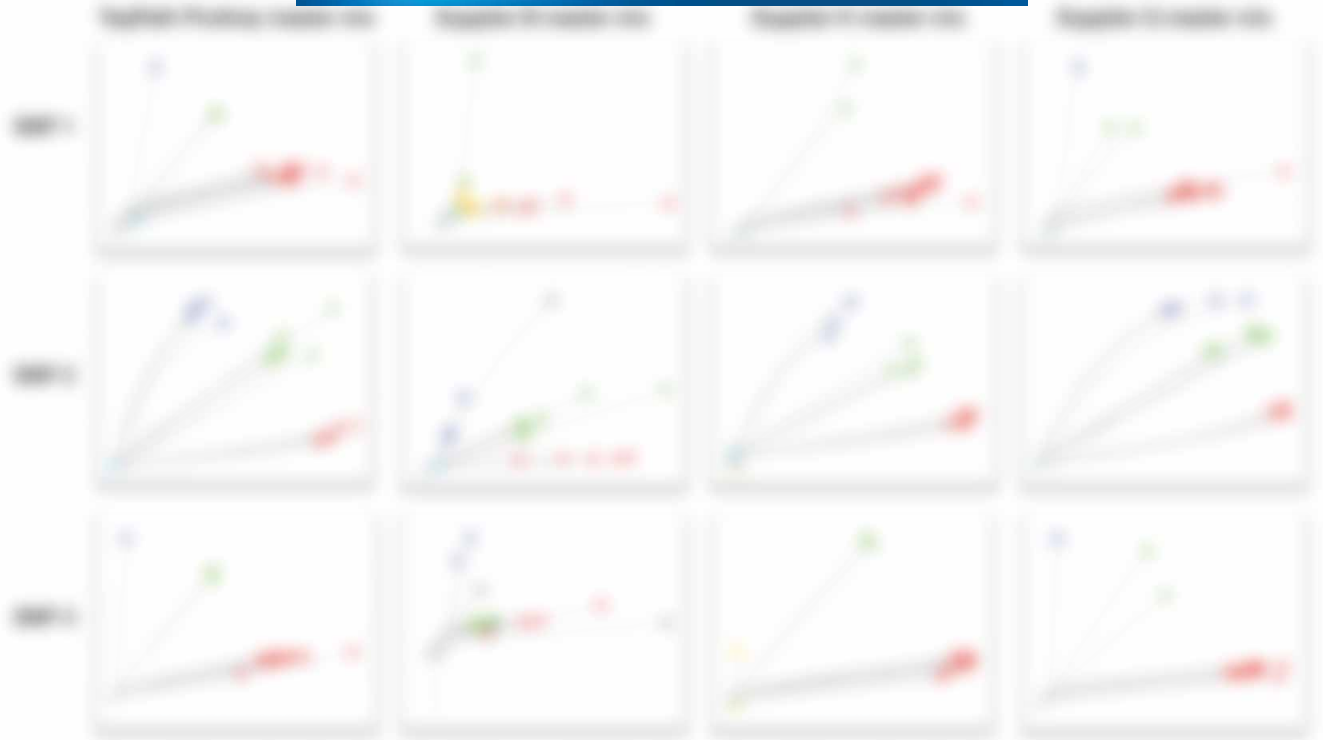
## Master mixes

The TaqPath ProAmp Multiplex Master Mix and other master mixes are designed to provide consistent and accurate results across multiple assays. The TaqPath ProAmp Multiplex Master Mix is a high-performance master mix that provides excellent cluster resolution and accurate genotype calls across multiple assays. The TaqPath ProAmp Multiplex Master Mix is designed to provide consistent and accurate results across multiple assays. The TaqPath ProAmp Multiplex Master Mix is a high-performance master mix that provides excellent cluster resolution and accurate genotype calls across multiple assays.

The TaqPath ProAmp Multiplex Master Mix is a high-performance master mix that provides excellent cluster resolution and accurate genotype calls across multiple assays. The TaqPath ProAmp Multiplex Master Mix is designed to provide consistent and accurate results across multiple assays. The TaqPath ProAmp Multiplex Master Mix is a high-performance master mix that provides excellent cluster resolution and accurate genotype calls across multiple assays.

**Thanks for reading!**  
Access the unabridged PDF and more at the QualTrak resources page.

[Keep reading](#)



**Figure 2. Genotyping results from TaqPath ProAmp Multiplex Master Mix and master mixes from other suppliers.** In three TaqMan™ genotyping assays performed under standard conditions, TaqPath ProAmp master mix consistently produced accurate genotype calls and excellent cluster resolution across multiple assays, demonstrating higher data precision compared to other suppliers' mixes.

## Predesigned and custom assay formats

ThermoFisher offers a wide range of assay formats to meet your needs. Our experts can help you design and optimize your assays for maximum efficiency and accuracy. We offer a variety of assay formats, including real-time PCR, digital PCR, and microfluidics. Our experts can help you choose the right assay format for your application and provide you with the necessary reagents and protocols.

For a complete list of assay formats and protocols, visit our website at [thermo.com/qpcr/biopharma](#). You can also contact our technical support team for more information.

ThermoFisher's assay formats are designed to be easy to use and highly reproducible. Our experts can help you troubleshoot any issues you may encounter and ensure that your results are accurate and reliable.

ThermoFisher's assay formats are also highly flexible, allowing you to adapt them to your specific needs. Our experts can help you modify your assays to accommodate different sample types and concentrations.

ThermoFisher's assay formats are also highly scalable, allowing you to run large numbers of samples. Our experts can help you optimize your assays for high-throughput screening and ensure that your results are consistent across all samples.

ThermoFisher's assay formats are also highly sensitive, allowing you to detect even low levels of target. Our experts can help you optimize your assays for maximum sensitivity and ensure that your results are accurate and reliable.

ThermoFisher's assay formats are also highly accurate, allowing you to obtain precise results. Our experts can help you optimize your assays for maximum accuracy and ensure that your results are consistent across all samples.

ThermoFisher's assay formats are also highly reproducible, allowing you to obtain consistent results. Our experts can help you optimize your assays for maximum reproducibility and ensure that your results are accurate and reliable.

**Thanks for reading!**  
Access the unabridged PDF and more at the QualTrak resources page.

[Keep reading](#)



## Ensure quality assurance (QA), control, and compliance with advanced technology and compliant software, documentation, and facilities

ThermoFisher offers a wide range of advanced technology and compliant software, documentation, and facilities to ensure quality assurance, control, and compliance. Our experts can help you choose the right technology and software for your application and provide you with the necessary reagents and protocols.

ThermoFisher's advanced technology and compliant software, documentation, and facilities are designed to be easy to use and highly reproducible. Our experts can help you troubleshoot any issues you may encounter and ensure that your results are accurate and reliable.

ThermoFisher's advanced technology and compliant software, documentation, and facilities are also highly flexible, allowing you to adapt them to your specific needs. Our experts can help you modify your assays to accommodate different sample types and concentrations.

## Leverage a robust supply chain to ensure uninterrupted PCR operations

ThermoFisher offers a wide range of advanced technology and compliant software, documentation, and facilities to ensure quality assurance, control, and compliance. Our experts can help you choose the right technology and software for your application and provide you with the necessary reagents and protocols.

ThermoFisher's advanced technology and compliant software, documentation, and facilities are designed to be easy to use and highly reproducible. Our experts can help you troubleshoot any issues you may encounter and ensure that your results are accurate and reliable.

ThermoFisher's advanced technology and compliant software, documentation, and facilities are also highly flexible, allowing you to adapt them to your specific needs. Our experts can help you modify your assays to accommodate different sample types and concentrations.

## Conclusions

Choosing to adopt better PCR technology and the right partner offers tremendous benefits for discovery, development, and manufacturing of biologics, including vaccines, cell and gene therapies, mAbs, and their biosimilars. To unlock the most effective instruments, master mixes, assays, and compliant software for your biologics program, find a vendor with a complete and thriving qPCR ecosystem.

Thermo Fisher is a long-established, trusted leader that can provide the leading cGMP-compliant QualTrak Real-Time PCR technologies and support to guide you through every PCR step of biologics development. Thermo Fisher guarantees continuous automation, accuracy, consistency, and compliance to its partners—setting them up for success from the first phase of discovery through widespread distribution of their biologic.

Visit [thermofisher.com/qpcr/biopharma](https://thermofisher.com/qpcr/biopharma) to get started with Thermo Fisher Scientific as your full-service biopharmaceutical PCR products provider. Our Applied Biosystems QualTrak biopharma-specific qPCR and dPCR products and workflows are designed to fast-track your biologics development pipeline, allowing you to develop the highest quality mAbs, biosimilars, vaccines, and other therapeutics in the shortest amount of time.

## References

1. Lu RM, Hwang YC, Liu IJ, et. al. (2020). Development of therapeutic antibodies for the treatment of diseases. *Journal of Biomedical Science* 27(1).
2. Chen Y, Monnard A, Santos da Silva J (7 June 2021). "[An inflection point for biosimilars.](#)" *McKinsey & Company*.
3. [Global Cell and Gene Therapy Market Report 2020: Market to recover and grow at a CAGR of 24.1% in 2023.](#) (7 January 2020) *Business Wire*.
4. Campbell H (26 August 2015). "[Video & infographic: Developing a new drug is actually harder than rocket science.](#)" *PhRMA*.

Learn more at [thermofisher.com/qpcr/biopharma](https://thermofisher.com/qpcr/biopharma)

**applied biosystems**