# Quick start guide: QuantStudio Absolute Q DNA Digital PCR Master Mix

This quick start guide provides concise instructions for preparing a digital PCR (dPCR) reaction for the Applied Biosystems<sup>™</sup> QuantStudio<sup>™</sup> Absolute Q<sup>™</sup> Digital PCR System. For detailed instructions on preparing and performing a dPCR experiment, refer to the QuantStudio<sup>™</sup> Absolute Q<sup>™</sup> Digital PCR Installation, Use, and Maintenance Guide (Pub. No. MAN0025621).

#### Getting started

- QuantStudio Absolute Q MAP16
  Digital PCR Plate
- 5 gaskets (per plate)
- Absolute Q DNA Digital PCR Master Mix (5X)
- QuantStudio Absolute Q
  Isolation Buffer
- Nuclease-free water
- Low-retention pipette tips
- Tabletop centrifuge
- Digital PCR assay

	Ft. I	V. I	V. I
Reagent	Final concentration	Volume per reaction	Volume per reaction with 10% overage*
Water	-	Fill to 9 μL	Fill to 10 μL
Absolute Q DNA Digital PCR Master Mix (5X)	1X	1.8 μL	2 μL
Digital PCR assay (40X or 20X)	1X	0.23 μL (40X) or 0.45 μL (20X)	0.25 μL (40X) or 0.50 μL (20X)
DNA sample	1–11,000 copies/μL**	Variable	Variable
Total	-	9 μL	10 μL

<sup>\*</sup> After calculating the number of reactions required, prepare dPCR mix for the appropriate number of reactions and scale those components by 10% for overage. Dilute assay accordingly to avoid pipetting less than 1  $\mu$ L volumes.

#### Prepare digital PCR reactions



Vortex the master mix well and combine the reagents shown in the table above in the order listed.



Mix dPCR reagents well.



Centrifuge at 10,000 x g for 1 minute (rotor centrifuge).

#### Load the QuantStudio Absolute Q MAP16 plate

Handle the MAP plate by its frame and work on a level, dust-free surface. Proceed to step 4, using a new pipette tip for each well.



At a 45° angle, load 9  $\mu L$  of the dPCR mixture to the bottom of the well. Pipette dPCR mixture only to the first stop.



At a 45° angle, load 15  $\mu L$  of Absolute Q Isolation Buffer. Pipette only to the first stop.



Apply 5 gaskets to the MAP plate. Orient the gaskets such that the "A" tail is on top of the plate.

### Setting up the run



Select the columns of the MAP plate that have been loaded.



Select "Start" to begin dPCR run



Define PCR thermal parameters. Set preheat step to 96°C for 10 minutes. No post-cycling steps required or recommended.



Select channel dye from the available options.

## Find out more at thermofisher.com/absoluteq



<sup>\*\*</sup> DNA copy number and dilution calculator can be found at thermofisher.com/dna-calculator