

QuantiGene Incubator Temperature Validation Kit

Cat No. QS0517

About QuantiGene Incubator Temperature Validation Kit

The QuantiGene® Incubator Temperature Validation Kit uses a National Institute of Standards and Technology (NIST) calibrated thermometer to assess the accuracy of the temperature of the oven incubator or shaking incubator used in all QuantiGene Assays. For the best assay results, we highly recommend verifying the incubator using the QuantiGene Incubator Temperature Validation Kit.

Intended Use

The QuantiGene Incubation Temperature Validation Kit is designed to precisely measure the temperatures in the incubators used in all QuantiGene Singleplex and QuantiGene Plex Assays.

Contents and Storage

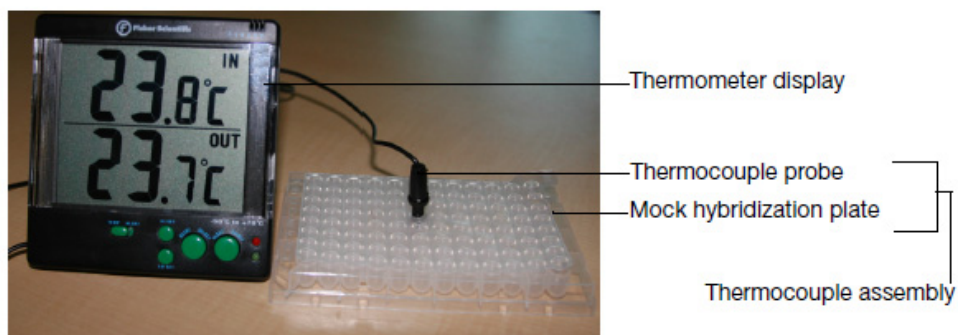
Component	Description	Quantity	Storage
Mock Hybridization Plate	A sealed microplate with a hole over well D7	1	15–30 °C
Thermometer	NIST calibrated digital thermometer with thermocouple probe	1	15–30 °C

Assembly Instructions

1. Activate the battery by removing the battery cover from the back of the thermometer. Two temperature readouts will appear on the thermometer display:
Out - Displays the current temperature of the thermocouple probe.
In - Displays the current temperature of the display environment.

NOTE: The date for recalibration is listed on the back of each thermometer.

2. Insert the thermometer probe into well D7 of the mock hybridization plate, ensuring that the probe is touching the bottom of the well.



Measuring the Incubator Temperature

1. Place the thermocouple assembly inside the pre-warmed incubator where you will place your QuantiGene assay plate and the thermometer display outside of the incubator. Ideally, the thermocouple assembly should be allowed to sit in the incubator overnight to reach equilibrium. Leave the thermometer for at least 2 hours before recording the measurement.
2. Measure and record the temperature of the incubator in different locations to ensure the temperature uniformity within the incubator. Temperature can vary by several degrees at different locations inside some incubators.
3. If necessary, adjust the incubator temperature so that the thermometer displays the assay-specified temperature.
4. Once the oven has been adjusted, check the temperature of the incubator at least once a month.

NOTE: The thermocouple/mock hybridization plate assembly can be left in the incubator for continual monitoring of the incubator temperature.

Measuring the VorTemp Shaking Incubator Temperature

1. Place an inverted plate lid into each of the two plate carriers in the pre-warmed VorTemp shaking incubator.
NOTE: Always use an inverted plate lid in the VorTemp when performing QuantiGene Plex Assays.
2. Place the thermocouple assembly on top of the inverted plate lid in one of the plate carriers. Place an empty Filter Plate with a plate seal on top of the other inverted Plate Lid in the plate carrier for balance.
3. Close the VorTemp's lid. Make sure to leave slack in the wire inside the VorTemp chamber so that the thermocouple probe will not get damaged during shaking.
4. Set the temperature and shaking speed specified in the QuantiGene Plex assay procedure and incubate overnight to permit the thermocouple assembly to reach equilibrium. If overnight incubation is not possible, incubate for at least 4 hours to ensure an accurate temperature reading.
5. Adjust the VorTemp shaking incubator temperature as necessary to match the assay-specified temperature displayed on the thermometer.

NOTE: The VorTemp display and the actual temperature measured by the thermometer might differ by 4 °C or more.

www.affymetrix.com/contact Please visit our website for international distributor contact information.
For research use only. Not for use in diagnostic procedures.