

QuantiGene® Incubator Temperature Validation Kit

Product Insert

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

Panomics' QuantiGene Incubation Temperature Validation Kit is designed to precisely measure the temperatures in the incubators used in all QuantiGene 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

Step	Action
1	<p>Activate the battery by removing the battery cover from the back of the thermometer. Two temperature readouts will appear on the thermometer display:</p> <p>Out—Displays the current temperature of the thermocouple probe.</p> <p>In—Displays the current temperature of the display environment.</p> <p>Note The date for recalibration is listed on the back of each thermometer.</p>
2	<p>Insert the thermometer probe into well D7 of the mock hybridization plate, ensuring that the probe is touching the bottom of the well.</p> <div style="text-align: right;">  <p>Thermometer display</p> <p>Thermocouple probe</p> <p>Mock hybridization plate</p> <p>Thermocouple assembly</p> </div>

Measuring the Incubator Temperature

Step	Action
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

Step	Action
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.

Contacting Panomics

For technical questions, please contact our technical support group by telephone at 1-877-726-6642 option 3 or email at techsupport@panomics.com (US and Canada). In Europe, contact techsupport_europe@panomics.com. In Asia Pacific, contact techsupport_asia@panomics.com. For an updated list of FAQs and product support literature, visit our website at www.panomics.com.

QuantiGene® is a registered trademark exclusively licensed to Panomics, Inc.

© 2008 Panomics, Inc. All rights reserved.

Panomics Inc., 6519 Dumbarton Circle, Fremont, CA 94555, Tel. (877) 726-6642