

QuantStudio™ 6 and 7 Flex Real-Time PCR Systems

Pub. No. MAN0019267 Rev. A.0

The *QuantStudio™ 6 and 7 Flex Real-Time PCR Systems Calibration and Performance Check Record* is a document designed to allow end users to perform and record the results of dye calibration and an instrument performance check. This document does not take the place of a qualification service performed by a certified Thermo Fisher Scientific field service engineer.

IMPORTANT! This document does not represent qualification of the QuantStudio™ 6 or 7 Flex Real-Time PCR System.

Customer and service information

Customer organization name:

Customer contact name:

Customer contact title:

Customer organization location:

Date performed:

Performed by:

Instrument information

Instrument model:

☐ QuantStudio™ 6 Flex ☐ QuantStudio™ 7 Flex

Instrument serial number:

Sample block and heated cover type:

☐ 96-Well ☐ Fast 96-Well ☐ 384-Well ☐ Array Card

Note: The Array Card block uses the 384-Well heated cover.

Sample block serial number:

Heated cover serial number:

Responsibilities

Customers should ensure that the *QuantStudio™ 6 and 7 Flex Real-Time PCR Systems Maintenance and Administration Guide* (Pub. No. 4489821 (v1.5 or earlier) or MAN0018828 (v1.6 or later)) is available, and that the appropriate consumables (listed in each section) have been obtained.

Optical calibrations

✓	Task																																																																						
<input type="checkbox"/>	Obtain the appropriate calibration plates listed below.																																																																						
	<table border="1"> <thead> <tr> <th></th> <th>Fast 96-Well</th> <th>96-Well</th> <th>384-Well</th> <th>Array Card</th> </tr> </thead> <tbody> <tr> <td>ROI and background kit</td> <td><input type="checkbox"/> 4432426</td> <td><input type="checkbox"/> 4432364</td> <td><input type="checkbox"/> 4432320</td> <td>N/A</td> </tr> <tr> <td>Normalization plates kit^[1]</td> <td><input type="checkbox"/> 4432432</td> <td><input type="checkbox"/> 4432370</td> <td><input type="checkbox"/> 4432308</td> <td>N/A</td> </tr> <tr> <td>FAM™ dye plate^[1]</td> <td><input type="checkbox"/> 4432389</td> <td><input type="checkbox"/> 4432327</td> <td><input type="checkbox"/> 4432271</td> <td>N/A</td> </tr> <tr> <td>TAMRA™ dye plate^[1]</td> <td><input type="checkbox"/> 4432414</td> <td><input type="checkbox"/> 4432352</td> <td><input type="checkbox"/> 4432296</td> <td>N/A</td> </tr> <tr> <td>ROX™ dye plate^[1]</td> <td><input type="checkbox"/> 4432402</td> <td><input type="checkbox"/> 4432340</td> <td><input type="checkbox"/> 4432284</td> <td>N/A</td> </tr> <tr> <td>VIC™ dye plate^[1]</td> <td><input type="checkbox"/> 4432396</td> <td><input type="checkbox"/> 4432334</td> <td><input type="checkbox"/> 4432278</td> <td>N/A</td> </tr> <tr> <td>NED™ dye plate^[1]</td> <td><input type="checkbox"/> 4432420</td> <td><input type="checkbox"/> 4432358</td> <td><input type="checkbox"/> 4432302</td> <td>N/A</td> </tr> <tr> <td>SYBR GREEN™ dye plate^[1]</td> <td><input type="checkbox"/> 4432408</td> <td><input type="checkbox"/> 4432346</td> <td><input type="checkbox"/> 4432290</td> <td>N/A</td> </tr> <tr> <td>ABY™ dye plate^[1]</td> <td><input type="checkbox"/> A24734</td> <td><input type="checkbox"/> A24738</td> <td><input type="checkbox"/> A24736</td> <td>N/A</td> </tr> <tr> <td>JUN™ dye plate^[1]</td> <td><input type="checkbox"/> A24735</td> <td><input type="checkbox"/> A24737</td> <td><input type="checkbox"/> A24733</td> <td>N/A</td> </tr> <tr> <td>MUSTANG PURPLE™ dye plate^[1]</td> <td><input type="checkbox"/> 4457328</td> <td><input type="checkbox"/> 4461599</td> <td><input type="checkbox"/> 4457334</td> <td>N/A</td> </tr> <tr> <td>10-dye spectral calibration kit^[2]</td> <td><input type="checkbox"/> A26342</td> <td><input type="checkbox"/> A26343</td> <td><input type="checkbox"/> A26341</td> <td>N/A</td> </tr> <tr> <td>Array Card calibration kit</td> <td>N/A</td> <td>N/A</td> <td>N/A</td> <td><input type="checkbox"/> 4432314</td> </tr> </tbody> </table>		Fast 96-Well	96-Well	384-Well	Array Card	ROI and background kit	<input type="checkbox"/> 4432426	<input type="checkbox"/> 4432364	<input type="checkbox"/> 4432320	N/A	Normalization plates kit ^[1]	<input type="checkbox"/> 4432432	<input type="checkbox"/> 4432370	<input type="checkbox"/> 4432308	N/A	FAM™ dye plate ^[1]	<input type="checkbox"/> 4432389	<input type="checkbox"/> 4432327	<input type="checkbox"/> 4432271	N/A	TAMRA™ dye plate ^[1]	<input type="checkbox"/> 4432414	<input type="checkbox"/> 4432352	<input type="checkbox"/> 4432296	N/A	ROX™ dye plate ^[1]	<input type="checkbox"/> 4432402	<input type="checkbox"/> 4432340	<input type="checkbox"/> 4432284	N/A	VIC™ dye plate ^[1]	<input type="checkbox"/> 4432396	<input type="checkbox"/> 4432334	<input type="checkbox"/> 4432278	N/A	NED™ dye plate ^[1]	<input type="checkbox"/> 4432420	<input type="checkbox"/> 4432358	<input type="checkbox"/> 4432302	N/A	SYBR GREEN™ dye plate ^[1]	<input type="checkbox"/> 4432408	<input type="checkbox"/> 4432346	<input type="checkbox"/> 4432290	N/A	ABY™ dye plate ^[1]	<input type="checkbox"/> A24734	<input type="checkbox"/> A24738	<input type="checkbox"/> A24736	N/A	JUN™ dye plate ^[1]	<input type="checkbox"/> A24735	<input type="checkbox"/> A24737	<input type="checkbox"/> A24733	N/A	MUSTANG PURPLE™ dye plate ^[1]	<input type="checkbox"/> 4457328	<input type="checkbox"/> 4461599	<input type="checkbox"/> 4457334	N/A	10-dye spectral calibration kit ^[2]	<input type="checkbox"/> A26342	<input type="checkbox"/> A26343	<input type="checkbox"/> A26341	N/A	Array Card calibration kit	N/A	N/A	N/A	<input type="checkbox"/> 4432314
	Fast 96-Well	96-Well	384-Well	Array Card																																																																			
ROI and background kit	<input type="checkbox"/> 4432426	<input type="checkbox"/> 4432364	<input type="checkbox"/> 4432320	N/A																																																																			
Normalization plates kit ^[1]	<input type="checkbox"/> 4432432	<input type="checkbox"/> 4432370	<input type="checkbox"/> 4432308	N/A																																																																			
FAM™ dye plate ^[1]	<input type="checkbox"/> 4432389	<input type="checkbox"/> 4432327	<input type="checkbox"/> 4432271	N/A																																																																			
TAMRA™ dye plate ^[1]	<input type="checkbox"/> 4432414	<input type="checkbox"/> 4432352	<input type="checkbox"/> 4432296	N/A																																																																			
ROX™ dye plate ^[1]	<input type="checkbox"/> 4432402	<input type="checkbox"/> 4432340	<input type="checkbox"/> 4432284	N/A																																																																			
VIC™ dye plate ^[1]	<input type="checkbox"/> 4432396	<input type="checkbox"/> 4432334	<input type="checkbox"/> 4432278	N/A																																																																			
NED™ dye plate ^[1]	<input type="checkbox"/> 4432420	<input type="checkbox"/> 4432358	<input type="checkbox"/> 4432302	N/A																																																																			
SYBR GREEN™ dye plate ^[1]	<input type="checkbox"/> 4432408	<input type="checkbox"/> 4432346	<input type="checkbox"/> 4432290	N/A																																																																			
ABY™ dye plate ^[1]	<input type="checkbox"/> A24734	<input type="checkbox"/> A24738	<input type="checkbox"/> A24736	N/A																																																																			
JUN™ dye plate ^[1]	<input type="checkbox"/> A24735	<input type="checkbox"/> A24737	<input type="checkbox"/> A24733	N/A																																																																			
MUSTANG PURPLE™ dye plate ^[1]	<input type="checkbox"/> 4457328	<input type="checkbox"/> 4461599	<input type="checkbox"/> 4457334	N/A																																																																			
10-dye spectral calibration kit ^[2]	<input type="checkbox"/> A26342	<input type="checkbox"/> A26343	<input type="checkbox"/> A26341	N/A																																																																			
Array Card calibration kit	N/A	N/A	N/A	<input type="checkbox"/> 4432314																																																																			
<input type="checkbox"/>	<p>Follow the instructions in the user guide to complete the ROI, background, and uniformity calibrations.</p> <p>ROI: <input type="checkbox"/> Passed</p> <p>Background: <input type="checkbox"/> Passed</p> <p>Uniformity: <input type="checkbox"/> Passed</p> <p>Note: If the background calibration fails, clean the block according to the instructions in the user guide.</p>																																																																						
<p>For Fast 96-Well, 96-Well, and 384-Well sample blocks with QuantStudio™ Software v1.5 or earlier only. If not applicable, select "N/A" below, then skip to the next section in this table.</p> <p><input type="checkbox"/> N/A</p>																																																																							
<input type="checkbox"/>	<p>Follow the instructions in the user guide to complete the desired pure dye calibrations.</p> <p>Note: To run the normalization calibration and the performance check, the instrument must be calibrated for FAM™, TAMRA™, ROX™, and VIC™ dyes at minimum.</p> <table border="0"> <thead> <tr> <th>Dye</th> <th>Result</th> <th>Dye</th> <th>Result</th> </tr> </thead> <tbody> <tr> <td>FAM™ dye</td> <td><input type="radio"/> Passed <input type="radio"/> Not calibrated</td> <td>SYBR GREEN™ dye</td> <td><input type="radio"/> Passed <input type="radio"/> Not calibrated</td> </tr> <tr> <td>TAMRA™ dye</td> <td><input type="radio"/> Passed <input type="radio"/> Not calibrated</td> <td>ABY™ dye</td> <td><input type="radio"/> Passed <input type="radio"/> Not calibrated</td> </tr> <tr> <td>ROX™ dye</td> <td><input type="radio"/> Passed <input type="radio"/> Not calibrated</td> <td>JUN™ dye</td> <td><input type="radio"/> Passed <input type="radio"/> Not calibrated</td> </tr> <tr> <td>VIC™ dye</td> <td><input type="radio"/> Passed <input type="radio"/> Not calibrated</td> <td>MUSTANG PURPLE™ dye</td> <td><input type="radio"/> Passed <input type="radio"/> Not calibrated</td> </tr> <tr> <td>NED™ dye</td> <td><input type="radio"/> Passed <input type="radio"/> Not calibrated</td> <td></td> <td></td> </tr> </tbody> </table>	Dye	Result	Dye	Result	FAM™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	SYBR GREEN™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	TAMRA™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	ABY™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	ROX™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	JUN™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	VIC™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	MUSTANG PURPLE™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	NED™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated																																																
Dye	Result	Dye	Result																																																																				
FAM™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	SYBR GREEN™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated																																																																				
TAMRA™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	ABY™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated																																																																				
ROX™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	JUN™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated																																																																				
VIC™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated	MUSTANG PURPLE™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated																																																																				
NED™ dye	<input type="radio"/> Passed <input type="radio"/> Not calibrated																																																																						
<input type="checkbox"/>	<p>Follow the instructions in the user guide to complete the normalization calibrations.</p> <p>FAM™ / ROX™ <input type="checkbox"/> Passed</p> <p>VIC™ / ROX™ <input type="checkbox"/> Passed</p>																																																																						

^[1] Not applicable to QuantStudio™ Software v1.6 or later.

^[2] Not applicable to QuantStudio™ Software v1.5 or earlier.

✓	Task												
For Fast 96-Well and 96-Well sample blocks with QuantStudio™ Software v1.6 or later only. If not applicable, select "N/A" below, then skip to the next section in this table.													
<input type="checkbox"/> N/A													
<input type="checkbox"/>	Follow the instructions in the user guide to complete the pure dye calibrations. <table border="1"> <thead> <tr> <th>Plate</th> <th>Dyes</th> <th>Result</th> </tr> </thead> <tbody> <tr> <td>Plate 1</td> <td>FAM™, VIC™, ROX™, SYBR GREEN™</td> <td><input type="checkbox"/> Passed</td> </tr> <tr> <td>Plate 2</td> <td>ABY™, JUN™, MUSTANG PURPLE™</td> <td><input type="checkbox"/> Passed</td> </tr> <tr> <td>Plate 3</td> <td>NED™, TAMRA™, Cy5™</td> <td><input type="checkbox"/> Passed</td> </tr> </tbody> </table>	Plate	Dyes	Result	Plate 1	FAM™, VIC™, ROX™, SYBR GREEN™	<input type="checkbox"/> Passed	Plate 2	ABY™, JUN™, MUSTANG PURPLE™	<input type="checkbox"/> Passed	Plate 3	NED™, TAMRA™, Cy5™	<input type="checkbox"/> Passed
Plate	Dyes	Result											
Plate 1	FAM™, VIC™, ROX™, SYBR GREEN™	<input type="checkbox"/> Passed											
Plate 2	ABY™, JUN™, MUSTANG PURPLE™	<input type="checkbox"/> Passed											
Plate 3	NED™, TAMRA™, Cy5™	<input type="checkbox"/> Passed											
For 384-Well sample block with QuantStudio™ Software v1.6 or later only. If not applicable, select "N/A" below, then skip to the next section in this table.													
<input type="checkbox"/> N/A													
<input type="checkbox"/>	Follow the instructions in the user guide to complete the pure dye calibrations. <table border="1"> <thead> <tr> <th>Plate</th> <th>Dyes</th> <th>Result</th> </tr> </thead> <tbody> <tr> <td>Plate 1</td> <td>FAM™, VIC™, ROX™, TAMRA™, SYBR GREEN™</td> <td><input type="checkbox"/> Passed</td> </tr> <tr> <td>Plate 2</td> <td>ABY™, JUN™, MUSTANG PURPLE™, NED™, Cy5™</td> <td><input type="checkbox"/> Passed</td> </tr> </tbody> </table>	Plate	Dyes	Result	Plate 1	FAM™, VIC™, ROX™, TAMRA™, SYBR GREEN™	<input type="checkbox"/> Passed	Plate 2	ABY™, JUN™, MUSTANG PURPLE™, NED™, Cy5™	<input type="checkbox"/> Passed			
Plate	Dyes	Result											
Plate 1	FAM™, VIC™, ROX™, TAMRA™, SYBR GREEN™	<input type="checkbox"/> Passed											
Plate 2	ABY™, JUN™, MUSTANG PURPLE™, NED™, Cy5™	<input type="checkbox"/> Passed											
For Array Card sample block only. If not applicable, select "N/A" below, then skip to the next section in this document.													
<input type="checkbox"/> N/A													
<input type="checkbox"/>	Follow the instructions in the user guide to complete the pure dye calibrations. <table border="1"> <tbody> <tr> <td>FAM™</td> <td><input type="checkbox"/> Passed</td> </tr> <tr> <td>ROX™</td> <td><input type="checkbox"/> Passed</td> </tr> <tr> <td>VIC™</td> <td><input type="checkbox"/> Passed</td> </tr> </tbody> </table>	FAM™	<input type="checkbox"/> Passed	ROX™	<input type="checkbox"/> Passed	VIC™	<input type="checkbox"/> Passed						
FAM™	<input type="checkbox"/> Passed												
ROX™	<input type="checkbox"/> Passed												
VIC™	<input type="checkbox"/> Passed												
<input type="checkbox"/>	Follow the instructions in the user guide to complete the normalization calibrations. <table border="1"> <tbody> <tr> <td>FAM™ / ROX™</td> <td><input type="checkbox"/> Passed</td> </tr> <tr> <td>VIC™ / ROX™</td> <td><input type="checkbox"/> Passed</td> </tr> </tbody> </table>	FAM™ / ROX™	<input type="checkbox"/> Passed	VIC™ / ROX™	<input type="checkbox"/> Passed								
FAM™ / ROX™	<input type="checkbox"/> Passed												
VIC™ / ROX™	<input type="checkbox"/> Passed												

Comments

If you have no comments, enter "Not applicable" below.

Performance check

✓	Task																																				
<input type="checkbox"/>	Obtain the appropriate RNase P plate listed below. <table border="1" style="width: 100%; border-collapse: collapse; margin-top: 10px;"> <thead> <tr> <th style="width: 35%;"></th> <th style="width: 15%;">Fast 96-Well</th> <th style="width: 15%;">96-Well</th> <th style="width: 15%;">384-Well</th> <th style="width: 20%;">Array Card</th> </tr> </thead> <tbody> <tr> <td>RNase P Instrument Verification Plate/Kit</td> <td style="text-align: center;"><input type="checkbox"/> 4351979</td> <td style="text-align: center;"><input type="checkbox"/> 4432382</td> <td style="text-align: center;"><input type="checkbox"/> 4455280</td> <td style="text-align: center;"><input type="checkbox"/> 4432265</td> </tr> </tbody> </table>		Fast 96-Well	96-Well	384-Well	Array Card	RNase P Instrument Verification Plate/Kit	<input type="checkbox"/> 4351979	<input type="checkbox"/> 4432382	<input type="checkbox"/> 4455280	<input type="checkbox"/> 4432265																										
	Fast 96-Well	96-Well	384-Well	Array Card																																	
RNase P Instrument Verification Plate/Kit	<input type="checkbox"/> 4351979	<input type="checkbox"/> 4432382	<input type="checkbox"/> 4455280	<input type="checkbox"/> 4432265																																	
For Fast 96-Well, 96-Well, and 384-Well sample blocks only. If not applicable, select "N/A" below, then skip to the next section in this table.																																					
<input type="checkbox"/>	N/A																																				
<input type="checkbox"/>	<p>Follow the instructions in the user guide to run the RNase P plate to complete the performance check. The instrument passes the installation specification if the inequality holds and the instrument successfully distinguishes between unknown populations A and B with a statistical confidence level of 99.7%. Performance is verified by the analysis equation below:</p> <p>$[(\text{CopyUnk}_2) - 3(\sigma_{\text{CopyUnk}_2})] > [(\text{CopyUnk}_1) + 3(\sigma_{\text{CopyUnk}_1})]$ where:</p> <ul style="list-style-type: none"> CopyUnk₂ = Average copy number (Qty Mean) of unknown population B (10000-copy population) $\sigma_{\text{CopyUnk}_2}$ = Standard deviation (Qty SD) of unknown population B (10000-copy population) CopyUnk₁ = Average copy number (Qty Mean) of unknown population A (5000-copy population) $\sigma_{\text{CopyUnk}_1}$ = Standard deviation (Qty SD) of unknown population A (5000-copy population) <p>Maximum number of outliers that can be removed:</p> <table border="1" style="width: 100%; border-collapse: collapse; margin-top: 10px;"> <thead> <tr> <th style="width: 35%;">Sample block</th> <th style="width: 15%;">Unknown outliers</th> <th style="width: 15%;">Standard outliers</th> <th style="width: 35%;">NTC outliers</th> </tr> </thead> <tbody> <tr> <td>Fast 96-Well or 96-Well</td> <td style="text-align: center;">6 per population</td> <td style="text-align: center;">1 per population</td> <td style="text-align: center;">0</td> </tr> <tr> <td>384-Well</td> <td style="text-align: center;">10 per population</td> <td style="text-align: center;">2 per population</td> <td style="text-align: center;">0</td> </tr> </tbody> </table> <p>When the run is complete, perform a primary analysis of the data and enter the values:</p> <table border="1" style="width: 100%; border-collapse: collapse; margin-top: 10px;"> <tbody> <tr> <td style="width: 33%;">CopyUnk₂ (Qty Mean 10K)</td> <td style="width: 33%;">$\sigma_{\text{CopyUnk}_2}$ (Qty SD 10K)</td> <td style="width: 34%;">Unknown B (10K) outliers</td> </tr> <tr> <td><input style="width: 90%;" type="text"/></td> <td><input style="width: 90%;" type="text"/></td> <td><input style="width: 90%;" type="text"/></td> </tr> <tr> <td>CopyUnk₁ (Qty Mean 5K)</td> <td>$\sigma_{\text{CopyUnk}_1}$ (Qty SD 5K)</td> <td>Unknown A (5K) outliers</td> </tr> <tr> <td><input style="width: 90%;" type="text"/></td> <td><input style="width: 90%;" type="text"/></td> <td><input style="width: 90%;" type="text"/></td> </tr> <tr> <td>Standard (1.25K) outliers</td> <td>Standard (2.5K) outliers</td> <td>Standard (5K) outliers</td> </tr> <tr> <td><input style="width: 90%;" type="text"/></td> <td><input style="width: 90%;" type="text"/></td> <td><input style="width: 90%;" type="text"/></td> </tr> <tr> <td>Standard (10K) outliers</td> <td>Standard (20K) outliers</td> <td>NTC outliers</td> </tr> <tr> <td><input style="width: 90%;" type="text"/></td> <td><input style="width: 90%;" type="text"/></td> <td><input style="width: 90%;" type="text"/></td> </tr> </tbody> </table> <p>Performance check: <input type="checkbox"/> Passed</p>	Sample block	Unknown outliers	Standard outliers	NTC outliers	Fast 96-Well or 96-Well	6 per population	1 per population	0	384-Well	10 per population	2 per population	0	CopyUnk ₂ (Qty Mean 10K)	$\sigma_{\text{CopyUnk}_2}$ (Qty SD 10K)	Unknown B (10K) outliers	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	CopyUnk ₁ (Qty Mean 5K)	$\sigma_{\text{CopyUnk}_1}$ (Qty SD 5K)	Unknown A (5K) outliers	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	Standard (1.25K) outliers	Standard (2.5K) outliers	Standard (5K) outliers	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	Standard (10K) outliers	Standard (20K) outliers	NTC outliers	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>
Sample block	Unknown outliers	Standard outliers	NTC outliers																																		
Fast 96-Well or 96-Well	6 per population	1 per population	0																																		
384-Well	10 per population	2 per population	0																																		
CopyUnk ₂ (Qty Mean 10K)	$\sigma_{\text{CopyUnk}_2}$ (Qty SD 10K)	Unknown B (10K) outliers																																			
<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>																																			
CopyUnk ₁ (Qty Mean 5K)	$\sigma_{\text{CopyUnk}_1}$ (Qty SD 5K)	Unknown A (5K) outliers																																			
<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>																																			
Standard (1.25K) outliers	Standard (2.5K) outliers	Standard (5K) outliers																																			
<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>																																			
Standard (10K) outliers	Standard (20K) outliers	NTC outliers																																			
<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>	<input style="width: 90%;" type="text"/>																																			

✓	Task																																
For Array Card sample block only. If not applicable, select "N/A" below, then skip to the next section in this document.																																	
<input type="checkbox"/> N/A																																	
<input type="checkbox"/>	<p>Follow the instructions in the user guide to prepare and run the RNase P kit to complete the performance check. The instrument passes the installation specification if the inequality holds and the instrument successfully distinguishes between unknown populations A and B with a statistical confidence level of 99.7%. Performance is verified by the analysis equation below:</p> <p>$[(\text{CopyUnk}_2) - 3(\sigma_{\text{CopyUnk}_2})] > [(\text{CopyUnk}_1) + 3(\sigma_{\text{CopyUnk}_1})]$ where:</p> <ul style="list-style-type: none"> CopyUnk₂ = Average copy number (Qty Mean) of unknown population B (1600-copy population) $\sigma_{\text{CopyUnk}_2}$ = Standard deviation (Qty SD) of unknown population B (1600-copy population) CopyUnk₁ = Average copy number (Qty Mean) of unknown population A (800-copy population) $\sigma_{\text{CopyUnk}_1}$ = Standard deviation (Qty SD) of unknown population A (800-copy population) <p>Maximum number of outliers that can be removed:</p> <table border="1"> <thead> <tr> <th>Sample block</th> <th>Unknown outliers</th> <th>Standard outliers</th> <th>NTC outliers</th> </tr> </thead> <tbody> <tr> <td>Array Card</td> <td>4 per population</td> <td>4 per population</td> <td>0</td> </tr> </tbody> </table> <p>When the run is complete, perform a primary analysis of the data and enter the values:</p> <table border="1"> <tbody> <tr> <td>CopyUnk₂ (Qty Mean 1.6K)</td> <td>$\sigma_{\text{CopyUnk}_2}$ (Qty SD 1.6K)</td> <td>Unknown B (1.6K) outliers</td> </tr> <tr> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> </tr> <tr> <td>CopyUnk₁ (Qty Mean 800)</td> <td>$\sigma_{\text{CopyUnk}_1}$ (Qty SD 800)</td> <td>Unknown A (800) outliers</td> </tr> <tr> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> </tr> <tr> <td>Standard (200) outliers</td> <td>Standard (400) outliers</td> <td>Standard (800) outliers</td> </tr> <tr> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> </tr> <tr> <td>Standard (1.6K) outliers</td> <td>Standard (3.2K) outliers</td> <td>NTC outliers</td> </tr> <tr> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> </tr> </tbody> </table> <p>Performance check: <input type="checkbox"/> Passed</p>	Sample block	Unknown outliers	Standard outliers	NTC outliers	Array Card	4 per population	4 per population	0	CopyUnk ₂ (Qty Mean 1.6K)	$\sigma_{\text{CopyUnk}_2}$ (Qty SD 1.6K)	Unknown B (1.6K) outliers	<input type="text"/>	<input type="text"/>	<input type="text"/>	CopyUnk ₁ (Qty Mean 800)	$\sigma_{\text{CopyUnk}_1}$ (Qty SD 800)	Unknown A (800) outliers	<input type="text"/>	<input type="text"/>	<input type="text"/>	Standard (200) outliers	Standard (400) outliers	Standard (800) outliers	<input type="text"/>	<input type="text"/>	<input type="text"/>	Standard (1.6K) outliers	Standard (3.2K) outliers	NTC outliers	<input type="text"/>	<input type="text"/>	<input type="text"/>
Sample block	Unknown outliers	Standard outliers	NTC outliers																														
Array Card	4 per population	4 per population	0																														
CopyUnk ₂ (Qty Mean 1.6K)	$\sigma_{\text{CopyUnk}_2}$ (Qty SD 1.6K)	Unknown B (1.6K) outliers																															
<input type="text"/>	<input type="text"/>	<input type="text"/>																															
CopyUnk ₁ (Qty Mean 800)	$\sigma_{\text{CopyUnk}_1}$ (Qty SD 800)	Unknown A (800) outliers																															
<input type="text"/>	<input type="text"/>	<input type="text"/>																															
Standard (200) outliers	Standard (400) outliers	Standard (800) outliers																															
<input type="text"/>	<input type="text"/>	<input type="text"/>																															
Standard (1.6K) outliers	Standard (3.2K) outliers	NTC outliers																															
<input type="text"/>	<input type="text"/>	<input type="text"/>																															

Comments


If you have no comments, enter "Not applicable" below.

Approval

Customer Name:

Customer Signature:

Date:

 Life Technologies Holdings Pte Ltd | Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256
For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0019267

Revision	Date	Description
A.0	15 April 2020	New document.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Cy5 is a trademark of GE Healthcare.