

MSD[®] Performance Qualification Kits

S 600/600MM and SQ 120/120MM Performance Qualification Kit

R31QQ-3



www.mesoscale.com[®]

Meso Scale Discovery

S 600/600MM and SQ 120/120MM Performance Qualification Kit

For use with the following MSD instruments: MESO® SECTOR S 600MM, MESO SECTOR® S 600, MESO QuickPlex® SQ 120MM, and MESO QuickPlex SQ 120.

This package insert must be read in its entirety before using this product.

FOR RESEARCH USE ONLY.

NOT FOR USE IN DIAGNOSTIC PROCEDURES.

Meso Scale Discovery

A division of Meso Scale Diagnostics, LLC.

1601 Research Boulevard

Rockville, MD 20850-3173 USA

www.mesoscale.com

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, mesoscale.com, www.mesoscale.com, methodicalmind.com, www.methodicalmind.com, Booster Pack, DISCOVERY WORKBENCH, InstrumentLink, MESO, MesoSphere, Methodical Mind, MSD GOLD, MULTI-ARRAY, MULTI-SPOT, QuickPlex, ProductLink, SECTOR, SECTOR HTS, SECTOR PR, SULFO-TAG, TeamLink, TrueSensitivity, TURBO-BOOST, TURBO-TAG, N-PLEX, R-PLEX, S-PLEX, T-PLEX, U-PLEX, V-PLEX, MSD (design), MSD (luminous design), Methodical Mind (head logo), 96 WELL SMALL-SPOT (design), 96 WELL 1-, 4-, 7-, 9-, & 10-SPOT (designs), 384 WELL 1- & 4-SPOT (designs), N-PLEX (design), R-PLEX (design), S-PLEX (design), T-PLEX (design), U-PLEX (design), V-PLEX (design), It's All About U, Spot the Difference, The Biomarker Company, and The Methodical Mind Experience are trademarks and/or service marks owned by or licensed to Meso Scale Diagnostics, LLC. All other trademarks and service marks are the property of their respective owners.

© 2014, 2019, 2024 Meso Scale Diagnostics, LLC. All rights reserved.

Table of Contents

Introduction	4
Reagents Supplied	4
Additional Materials and Equipment	5
Safety	5
Test Protocol	6
Records and Results	6
Troubleshooting	8

Contact Information

Customer Service

Phone: 1-240-314-2795

Fax: 1-301-990-2776

Email: CustomerService@mesoscale.com

Scientific Support

Phone: 1-240-314-2798

Fax: 1-240-632-2219 attn: Scientific Support

Email: ScientificSupport@mesoscale.com

Introduction

This Performance Qualification (PQ) Kit provides a rapid and convenient method for verifying instrument performance. The qualification procedure measures consistency of signals across a plate and dark (electronic) noise of the instrument.

The entire PQ test should take approximately 25 to 30 minutes and should be performed using the MSD reagents provided in the PQ Kit. Only those components from the lots specified in the kit's certificate of analysis (COA) should be used for each test. Mixing components from different manufacturing lots may compromise performance.

Mean signal, standard deviation, and %CV values are calculated and compared with specifications. Dark noise is measured as a standard deviation of the signal values from wells in which no electrochemiluminescence is occurring (i.e. empty wells). The source of these signals is electronic noise in the analog to digital conversion.

All performance information, including standard signal levels, is based on 150 μ L read volume in 96-well MSD assay plates using standard read parameters in the specified operating environment.

The intervals for conducting the PQ are defined by the operator and the quality system of the lab in which the instrument is being used.

The PQ procedure can only verify that the instrument is performing within specifications.

Reagents Supplied

Product Description	Storage	Quantity per Kit
96-Well 1 Spot SECTOR™ Plate	Room Temperature	10 plates
Free Tag ECL 15,000 (250 mL)	Room Temperature	1 bottle
PQ Low Control (200 mL)	Room Temperature	1 bottle

A lot-specific COA is included in each kit.

Important:

- ☐ Testing should be performed at 20–26 °C.
- ☐ Recap reagent bottles immediately after use. Leaving bottles open may cause spills or evaporation that will alter the concentration of these reagents and may result in failure of the PQ.
- ☐ Extended exposure of MSD assay plates to the air can affect the measured signal. Immediately reseal unused plates in original packaging to prevent exposure to the air.
- ☐ Do not mix components from different kit lots.

Additional Materials and Equipment

- ❑ MSD electronic test plate (provided at the time of instrument installation).
- ❑ Multi-channel pipettor, capable of dispensing 150 µL/well into a 96-well plate (customer provided). Ensure that the pipettor is calibrated.

Safety

Consult the *MSD Reader Safety Guide* for safety precautions and regulations concerning the handling of materials and the instrument's electrical and mechanical components before working with MSD plate readers.

Free Tag ECL and PQ Low Control are chemical solutions and should be handled appropriately. Proper care should be taken to prevent spills. Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling kit components. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines.

Please review additional product-specific safety information found in the applicable safety data sheet(s) (SDS), which can be obtained from MSD Customer Service or at www.mesoscale.com.

Test Protocol

1. Confirm the following:
 - ☐ The testing environment (lab) is within the recommended operating temperature for the instrument: 20–26 °C.
 - ☐ The instrument is plugged into a power source and powered on.
 - ☐ The instrument computer is powered on and the operating system is fully initialized.
 - ☐ The DISCOVERY WORKBENCH® or Methodical Mind® software is running.
 - ☐ The instrument camera is at operating temperature.
2. Load the instrument's electronic test plate on the instrument's input stack and read the plate. Verify the operation of the instrument confirming the following wells have signals greater than 2,000 counts: A9, B10, C11, D12, H1, G2, F3, and E4, and all other wells have signals lower than 200 counts.

Note: Review the instrument user guide for instructions on how to log into the software and read an MSD plate.
3. Record the following items in the records table provided below:
 - ☐ Lot numbers and expiration dates of all reagents
 - ☐ Plate barcodes
 - ☐ The obtained mean signal for the Free Tag and the PQ Low Control from the PQ kit COA
 - ☐ Pipettor information
 - ☐ Lab temperature and comments, if any

Note: Confirm that the lot numbers match those listed on the COA for this PQ kit.
4. Prepare 3 plates as follows:
 - ☐ Fill plate #1 with 150 µL of Free Tag in each well.
 - ☐ Fill plate #2 with 150 µL of PQ Low Control in each well.
 - ☐ Leave plate #3 empty.

Note: Use reverse pipetting technique and pipette to the bottom corner of the well to avoid creating bubbles. Use a clean pipette tip to burst any bubbles.
5. Seal and incubate plates 1 and 2 for 15 minutes (\pm 30 seconds) at 20–26 °C. Record the incubation time in the records table.
6. Read plates 1, 2, and 3 on the instrument one at a time and record the appropriate details in the results table provided below.
 - ☐ Calculate for plate #1: Experimental mean signal and coefficient of variation (%CV) values for the Free Tag. Enter the values in the results table.
 - ☐ Calculate for plate #2: Experimental mean signal and standard deviation (SD) for the PQ Low Control. Enter the values in the results table.
 - ☐ Calculate for plate #3: Experimental mean signal and SD for the dark noise. Enter the values in the results table.
7. Compare the calculated results obtained in step 4 with the specifications in the results table.
 - ☐ If all results are within the acceptable range, the test result is PASS.
 - ☐ If any one of the results is outside of the acceptable range, refer to the Troubleshooting section for guidance.

Records and Results

Operator:_____

Date:_____

Instrument Serial Number:_____

RECORDS TABLE	Lot Number	Expiration Date	Obtained Mean Signal from the COA
Performance Qualification Kit			
Free Tag			
PQ Low Control			
	Barcode	Incubation Time (Minutes)	
Plate 1			
Plate 2			
Plate 3			
	ID/Serial Number	Calibration Due Date	
Multichannel Pipettor			
	Ambient Temperature	Within Range 20–26 °C (Y/N)? If no, please provide comments	
Lab Temperature			

RESULTS TABLE	Description	Experimental Mean Signal	Signal Specification		PASS/FAIL
			Minimum Value	Maximum Value	
Plate 1	Free Tag		12,000	18,000	
Plate 2	PQ Low Control		25	100	
Plate 3	Dark Noise		-16	16	
	Description	Calculated Value	%CV and SD Specification		PASS/FAIL
Plate 1	Free Tag %CV		≤ 6 %		
Plate 2	PQ Low Control SD		≤ 25 counts		
Plate 3	Dark Noise SD		≤ 16 counts		

Troubleshooting

If the performance qualification test does not meet specifications, repeat the test after confirming the following:

1. This PQ test should be performed in a lab where the temperature is within the operational range of the reader (20–26 °C). If the PQ kit was stored outside of this temperature range, reagents and plates should be equilibrated to 20–26 °C prior to use, as follows:
 - Incubate the test plates at 20–26 °C for ≥ 30 minutes.
 - Incubate Free Tag and PQ Low Control bottles in a 20–26 °C water bath for ≥ 30 minutes. Partially submerge the bottles in the water bath, keeping the cap of the bottles above water level.
2. Plates and reagents were properly stored.
 - Ensure Free Tag and PQ Low Control reagent bottles have been properly stored before use. Leaving bottles open may cause evaporation that will alter the concentration of these reagents and may result in failure of the PQ.
 - Extended exposure of MSD assay plates to the air can affect the measured signal. Ensure unused plates have been properly stored in original packaging.
3. The expiration dates of all plates and reagents are not past due.
4. The correct protocol was followed.
 - The incubation time of Free Tag should be 15 minutes (± 30 seconds).
 - The volume of reagent per well should be 150 μL .
 - There were no bubbles in the wells.
5. All lot numbers of plates and reagents used match those listed in the PQ kit COA.

If the specifications are not met when the PQ test is repeated, please contact ScientificSupport@mesoscale.com.