



## CERTIFICATE OF ANALYSIS

**Product Description:** S 600/600MM and SQ 120/120MM Performance Qualification Kit  
**Catalog Number:** R31QQ-3  
**Lot Number:** PQ300126  
**Manufacture Date:** 29 JUL 2019  
**Expiration Date:** 30 JUN 2021

**Statement:** This product was manufactured and tested according to approved MSD Standard Operating Procedures.

Components for Testing S 600/600MM and SQ 120/120MM Performance Qualification Kit	Catalog Number	Lot Number	Expiration Date
SECTOR Imager Plate	L15XA-3	21J2KA	30 JUN 2021
MSD Free Tag ECL 15000	N/A	Y0260163	28 FEB 2022
PQ Low Control	R92TE-2	Y0110074	31 MAY 2022

S 600/600MM and SQ 120/120MM Performance Qualification Kit Specifications			
Test	QC Obtained Mean Signal	Signal Specifications	
		Minimal Value	Maximal Value
MSD Free Tag ECL 15000	15,065	12,000	18,000
PQ Low Control	69	25	100
Dark Noise	2	-16	16
Test	%CV and SD Specification		
MSD Free Tag ECL 15000	0 % - 6 %		
PQ Low Control SD	0 - 25		
Dark Noise SD	0 - 16		

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality	<i>Karen Hamilla</i>	06 AUG 2019