



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S 600/600MM and SQ 120/120MM Performance Qualification Kit

Catalog Number: R31QQ-3

Lot Number: PQ300138

Expiration Date: 30 JUN 2022

Storage Temperature: Room Temperature

Components	Catalog Number	Lot Number	Expiration Date
SECTOR Imager Plate	L15XA	21K2BA	30 JUN 2022
MSD Free Tag ECL 15000	N/A	Y0260167	31 AUG 2023
PQ Low Control	R92TE	Y0110083	30 SEP 2023

Test	QC Obtained Mean Signal	Signal Specifications	
		Minimum Value	Maximum Value
MSD Free Tag ECL 15000	15416	12000	18000
PQ Low Control	79	25	100
Dark Noise	3	-16	16
Test	%CV and SD Specification		
MSD Free Tag ECL 15000	0-6.0%		
PQ Low Control SD	0-25		
Dark Noise SD	0-16		

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		02 OCT 2020