



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: PQ300137

Expiration Date: 31 MAY 2022

Storage Temperature: Room Temperature

Components	Catalog Number	Lot Number	Expiration Date
SECTOR Imager Plate	L15XA	21K1QA	31 MAY 2022
MSD Free Tag ECL 15000	N/A	Y0260166	31 MAY 2023
PQ Low Control	R92TE	Y0110081	30 APR 2023

Test	QC Obtained Mean Signal	Signal Specifications	
		Minimum Value	Maximum Value
MSD Free Tag ECL 15000	15177	12000	18000
PQ Low Control	70	25	100
Dark Noise	1	-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		15 JUN 2020