



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: PQ300134

Expiration Date: 31 JUL 2021

Storage Temperature: Room Temperature

Components	Catalog Number	Lot Number	Expiration Date
SECTOR Imager Plate	L15XA	21J35A	31 JUL 2021
MSD Free Tag ECL 15000	N/A	Y0260165	31 DEC 2022
PQ Low Control	R92TE	Y0110073	31 JAN 2022

Test	QC Obtained Mean Signal	Signal Specifications	
		Minimum Value	Maximum Value
MSD Free Tag ECL 15000	14653	12000	18000
PQ Low Control	72	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		14 FEB 2020