



### 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:** PQ300139

**Expiration Date:** 30 JUN 2021

**Storage Temperature:** Room Temperature

Components	Catalog Number	Lot Number	Expiration Date
SECTOR Imager Plate	L15XA	21J2DA	30 JUN 2021
MSD Free Tag ECL 15000	N/A	Y0260167	31 AUG 2023
PQ Low Control	R92TE	Y0110074	31 MAY 2022

Test	QC Obtained Mean Signal	Signal Specifications	
		Minimum Value	Maximum Value
MSD Free Tag ECL 15000	15056	12000	18000
PQ Low Control	62	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		02 SEP 2020