



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® SECTOR Imager Performance Qualification Kit

Catalog Number: R31QQ-3

Lot Number: PQ300130

Expiration Date: 31 OCT 2021

Storage Temperature: Room Temperature

Components	Catalog Number	Lot Number	Expiration Date
SECTOR Imager Plate	L15XA	21J4CA	31 OCT 2021
MSD Free Tag ECL 15000	N/A	Y0260164	30 SEP 2022
PQ Low Control	R92TE	Y0110074	31 MAY 2022

Test	QC Obtained Mean Signal	Signal Specifications	
		Minimum Value	Maximum Value
MSD Free Tag ECL 15000	14477	12 000	18 000
PQ Low Control	66	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		28 OCT 2019