



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® SECTOR Imager Performance Qualification Kit

Catalog #/Size: R31QQ-3

Lot Number: PQ300120

Expiration Date: 28 FEB 2021

Statement: This product was manufactured and tested according to approved MSD documents.

SECTOR Imager Plate					
Catalog Number	L15XA	Lot Number	21J0CA	Expiration Date	28 FEB 2021

Free Tag 15,000					
Catalog Number	N/A	Lot Number	Y0260160	Expiration Date	30 NOV 2021

PQ Low Control					
Catalog Number	R92TE	Lot Number	Y0110073	Expiration Date	31 JAN 2022

Specifications:	Max	Min
Free Tag Mean	18 000	12 000
Free Tag CV	6	0
PQ Low Control Mean	100	25
PQ Low Control SD	25	0
Dark Signal Mean	16	-16
Dark Noise (SD)	16	0

MSD QC Results:					
Using Instrument	Sector Imager		S/N	1201130115105	
Free TAG Mean	15 350	PQ Low Control Mean	79	Dark Signal Mean	2.9

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality		17 APR 2019