



CERTIFICATE OF ANALYSIS

Page 1 of 0

Product Description: MSD[®] SECTOR Imager Performance Qualification Kit

Catalog #/Size: R31QQ-3

Lot Number: PQ300088

Expiration Date: 31 MAY 2019

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

SECTOR Imager Plate					
Catalog Number	L15XA / L11XA	Lot Number	21H1VA	Expiration Date	31 MAY 2019

Free Tag 15,000					
Catalog Number	N/A	Lot Number	Y0260156	Expiration Date	31 MAY 2020

PQ Low Control					
Catalog Number	R92TE	Lot Number	Y0320010	Expiration Date	30 APR 2020

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	14 376	PQ Low Control Mean	66	Dark Signal Mean	2

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	20 JUN 2017