



### CERTIFICATE OF ANALYSIS

Page 2 of 1

**Product Description:** MSD® SECTOR Imager Performance Qualification Kit

**Catalog #/Size:** R31QQ-3

**Lot Number:** PQ300106

**Expiration Date:** 31 MAR 2020

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

SECTOR Imager Plate					
Catalog Number	L15XA / L11XA	Lot Number	21I0MA	Expiration Date	31 MAR 2020

Free Tag 15,000					
Catalog Number	N/A	Lot Number	Y0260158	Expiration Date	30 SEP 2020

PQ Low Control					
Catalog Number	R92TE	Lot Number	Y0320013	Expiration Date	28 FEB 2021

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	15 329	PQ Low Control Mean	73	Dark Signal Mean	5

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>	22 MAY 2018