



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® SECTOR Imager Performance Qualification Kit

Catalog #/Size: R31QQ-3

Lot Number: PQ300122

Expiration Date: 31 JAN 2021

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

SECTOR Imager Plate					
Catalog Number	L15XA	Lot Number	21J0IA	Expiration Date	31 JAN 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	15 519	PQ Low Control Mean	67	Dark Signal Mean	1

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