



## CERTIFICATE OF ANALYSIS

**Product Description:** MSD SECTOR® Imager Performance Qualification Kit  
**Catalog Number:** R31QQ-3  
**Lot Number:** PQ300127  
**Manufacture Date:** 22 JUL 2019  
**Expiration Date:** 30 JUN 2021

**Statement:** This product was manufactured and tested according to approved MSD Standard Operating Procedures.

Components for Testing MSD SECTOR® Imager Performance Qualification Kit	Catalog Number	Lot Number	Expiration Date
SECTOR Imager Plate	L15XA-3	21J2EA	30 JUN 2021
MSD Free Tag ECL 15000	N/A	Y0260163	28 FEB 2022
PQ Low Control	R92TE-2	Y0110074	31 MAY 2022

MSD SECTOR® Imager Performance Qualification Kit			
Test	QC Obtained Mean Signal	Signal Specifications	
		Minimal Value	Maximal Value
MSD Free Tag ECL 15000	15,582	12,000	18,000
PQ Low Control	69	25	100
Dark Noise	0.1	-16	16
Test	%CV and SD Specification		
MSD Free Tag ECL 15000	0 % - 6 %		
PQ Low Control SD	0 - 25		
Dark Noise SD	0 - 16		

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 NOV 2019