



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® SECTOR PR2 Model 1800 Performance Qualification Kit

Catalog #/Size: R61QQ-2

Lot Number: PQ400038

Expiration Date: 30 JUN 2020

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

SECTOR PR2 Plate					
Catalog Number	2072-0001	Lot Number	1VIOQA	Expiration Date	30 JUN 2020

Free Tag 15,000					
Catalog Number	N/A	Lot Number	Y0260159	Expiration Date	30 APR 2021

Specifications:	Max	Min
Free Tag Mean	20 000	10 000
Free Tag CV	15	0
Dark Signal Mean	16	-16
Dark Noise (SD)	40	0

MSD QC Results:			
Using Instrument	PR2	S/N	19000808200111
Free TAG Mean	14 136	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>	12 SEP 2018