



### CERTIFICATE OF ANALYSIS

Page 1 of 0

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

**Catalog #/Size:** R61QQ-2

**Lot Number:** PQ400037

**Expiration Date:** 30 APR 2020

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

SECTOR PR2 Plate					
<b>Catalog Number</b>	2072-0001	<b>Lot Number</b>	1VIOEA	<b>Expiration Date</b>	30 APR 2020

Free Tag 15,000					
<b>Catalog Number</b>	N/A	<b>Lot Number</b>	Y0260159	<b>Expiration Date</b>	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:			
<b>Using Instrument</b>	PR2	<b>S/N</b>	19000808200111
<b>Free TAG Mean</b>	14 293	<b>Dark Signal Mean</b>	-10

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	07 SEP 2018