



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

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

**Lot Number:** PQ400040

**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

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

MSD QC Results:			
<b>Using Instrument</b>	PR2	<b>S/N</b>	19000808200111
<b>Free TAG Mean</b>	13 663	<b>Dark Signal Mean</b>	-3

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

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	06 NOV 2018