



**RFE MEMO**

Page 1 of 1

**Product Description(s):** MSD® SECTOR PR Performance Qualification Kit


**Catalog Number(s) / Size(s):** R21QQ-2

**Lot Number:** PQ100027

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of PQ100027 has been extended to 30 Apr 2020.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
MULTI-ARRAY® PR Plate	01H01A	30 APR 2019	30 APR 2020

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar		20 MAR 2019

This memo supersedes all other statements regarding date of expiration for this product.



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® SECTOR PR Performance Qualification Kit

**Catalog #/Size:** R21QQ-2

**Lot Number:** PQ100027

**Expiration Date:** 30 APR 2019

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

SECTOR PR Plate					
<b>Catalog Number</b>	L13XA / L17XA	<b>Lot Number</b>	01H01A	<b>Expiration Date</b>	30 APR 2019

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

Read Buffer T(1X)					
<b>Catalog Number</b>	R92TA	<b>Lot Number</b>	Y0110067	<b>Expiration Date</b>	31 MAY 2020

Specifications:	Max	Min
Free Tag Mean	18 000	12 000
Free Tag CV	6	0
Read Buffer T(1X) Mean	300	25
Read Buffer T(1X) SD	40	0
Dark Signal Mean	16	-16
Dark Noise (SD)	40	0

MSD QC Results:					
<b>Using Instrument</b>	PR400		<b>S/N</b>	1150090325055	
<b>Free TAG Mean</b>	14 784	<b>Read Buffer T(1X) Mean</b>	61	<b>Dark Signal Mean</b>	1

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>	28 NOV 2018