



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® SECTOR PR Performance Qualification Kit

Catalog #/Size: R21QQ-2

Lot Number: PQ100026

Expiration Date: 28 FEB 2019

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

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

Free Tag 15,000					
Catalog Number	N/A	Lot Number	Y0260153	Expiration Date	28 FEB 2019

Read Buffer T(1X)					
Catalog Number	R92TA	Lot Number	Y0110066	Expiration Date	29 FEB 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:					
Using Instrument	PR400		S/N	1150090325055	
Free TAG Mean	15 164	Read Buffer T(1X) Mean	71	Dark Signal Mean	-1

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>	11 SEP 2017