



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® PR2 Model 1800 Performance Qualification Kit

Catalog #/Size: R61QQ-2

Lot Number: PQ400043

Expiration Date: 30 SEP 2020

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

SECTOR PR2 Plate					
Catalog Number	2072-0001	Lot Number	1VI0UA	Expiration Date	30 SEP 2020

Free Tag 15,000					
Catalog Number	N/A	Lot Number	Y0260160	Expiration Date	30 NOV 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	14 546	Dark Signal Mean	-2.6

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 FEB 2019