

CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog #/Size: R61QQ-2 Lot Number: PQ400036 Expiration Date: 31 JUL 2019

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

SECTOR PR2 Plate					
Catalog Number	2072-0001	Lot Number	1VH01A	Expiration Date	31 JUL 2019

Free Tag 15,000					
Catalog Number	N/A	Lot Number	Y0260158	Expiration Date	30 SEP 2020

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	16 540	Dark Signal Mean	-2	

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	Dawn Stollar	06 MAR 2018

COA-9181-A QA-FM-191-A