



Rev. B

CERTIFICATE OF ANALYSIS

COA – 04416r

Product Description: MSD[®] Performance Qualification Kit, Sector PR2, Model 1800**Catalog #/Size:** R61QQ-2**Lot Number:** PQ400022**Expiration Date:** 30 JUN 2016**Statement:** This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

SECTOR PR2 Plate Lot Number: 1VE0TA

SECTOR PR2 Plate Catalog Number: 3-2072-020001

SECTOR PR2 Plate Expiration Date: 30 JUN 2016

Free Tag 15,000 Lot Number: Y0260146

Free Tag 15,000 Catalog Number: N/A

Free Tag 15,000 Expiration Date: 31 JUL 2017

Specifications:

	Max	Min
Free Tag Mean	20 000	10 000
Free Tag CV %	15.0	0.00
Dark Signal Mean	16	-16
Dark Noise (SD)	40	0

MSD QC Results: Using instrument: PR2

S/N: 19000808200111

Free TAG Mean: 14 935

Dark Signal Mean: -4

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		19 SEP 2014
Review/Approval	Anastasia Belilovets	Quality Control		19 SEP 2014
Review/Approval	Richard Mann	Quality Assurance		19 SEP 2014