



CERTIFICATE OF ANALYSIS

COA – 05200

Product Description: MSD[®] Performance Qualification Kit, Sector PR2, Model 1800

Catalog #/Size: R61QQ-2

Lot Number: PQ400025

Expiration Date: 31 AUG 2016

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

SECTOR PR2 Plate Lot Number: 1VE1BA

SECTOR PR2 Plate Catalog Number: 3-2072-020001

SECTOR PR2 Plate Expiration Date: 31 AUG 2016

Free Tag 15,000 Lot Number: Y0260150

Free Tag 15,000 Catalog Number: N/A

Free Tag 15,000 Expiration Date: 31 MAY 2018

Specifications:

	Max	Min
Free Tag Mean	18 000	12 000
Free Tag CV %	10.0	0.00
Dark Signal Mean	16	-16
Dark Noise (SD)	40	0




MSD QC Results: Using instrument: PR2

S/N: 19000808200111

Free TAG Mean: 14 786

Dark Signal Mean: -2

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		16 JUL 2015
Review/Approval	Anastasia Belilovets	Quality Assurance		16 JUL 2015
Review/Approval	Michael Grossman	Quality Assurance		16 JUL 2015