



CERTIFICATE OF ANALYSIS

COA – 04349

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

Catalog #/Size: R61QQ-2

Lot Number: PQ400020

Expiration Date: 30 APR 2016

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

SECTOR PR2 Plate Lot Number: 1VE0PA

SECTOR PR2 Plate Catalog Number: 2072-0001

SECTOR PR2 Plate Expiration Date: 30 APR 2016

Free Tag 15,000 Lot Number: Y0260145

Free Tag 15,000 Catalog Number: N/A

Free Tag 15,000 Expiration Date: 31 JAN 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: 17 850

Dark Signal Mean: -6

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		12 JUN 2014
Review/Approval	Dana Short	Quality Control		12 JUN 2014
Review/Approval	Richard Mann	Quality Assurance		12 JUN 2014