



CERTIFICATE OF ANALYSIS

COA – 04689

Product Description: MSD[®] Performance Qualification Kit, Sector Imager

Catalog #/Size: R31QQ-3

Lot Number: PQ300046

Expiration Date: 31 OCT 2016

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

SECTOR Imager Plate Lot Number: 21E2TA

SECTOR Imager Plate Catalog Number: L15XA

SECTOR Imager Plate Expiration Date: 31 OCT 2016

Free Tag 15,000 Lot Number: Y0260147

Free Tag 15,000 Catalog Number: N/A

Free Tag 15,000 Expiration Date: 31 OCT 2017

PQ Low Control Lot Number: Y0320005

PQ Low Control Catalog Number: R92TE

PQ Low Control Expiration Date: 30 JUN 2017

Specifications:

	Max	Min
Free Tag Mean	18 000	12 000
Free Tag CV	6	0
PQ Low Control Mean	100	25
PQ Low Control Mean SD	25	0
Dark Signal Mean	16	-16
Dark Noise (SD)	16	0

MSD QC Results: Using instrument SI: 6000 S/N: 1201130115105

Free TAG Mean: 15 528

PQ Low Control Mean: 61

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		18 DEC 2014
Review/Approval	Anastasia Belilovets	Quality Control		18 DEC 2014
Review/Approval	Richard Mann	Quality Assurance		18 DEC 2014