



### CERTIFICATE OF ANALYSIS

COA – 04524

**Product Description:** MSD® Performance Qualification Kit, Sector Imager

**Catalog #/Size:** R31QQ-3

**Lot Number:** PQ300045

**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: 21E2QA  
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: 15 137  
PQ Low Control Mean: 67  
Dark Signal Mean: -2

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

	Name	Function	Signature	Date
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		12 NOV 2014
<b>Review/Approval</b>	Anastasia Belilovets	Quality Control		12 NOV 2014
<b>Review/Approval</b>	Richard Mann	Quality Assurance		12Nov2014