



CERTIFICATE OF ANALYSIS

COA – 04027

Product Description: MSD® Performance Qualification Kit, Sector Imager

Catalog #/Size: R31QQ-3

Lot Number: PQ300034

Expiration Date: 30 JUN 2015

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

SECTOR Imager Plate Lot Number: 21D2CA
SECTOR Imager Plate Catalog Number: L15XA
SECTOR Imager Plate Expiration Date: 30 JUN 2015
Free Tag 15,000 Lot Number: Y0260144
Free Tag 15,000 Catalog Number: N/A
Free Tag 15,000 Expiration Date: 30 NOV 2016
PQ Low Control Lot Number: Y0320001
PQ Low Control Catalog Number: R92TE
PQ Low Control Expiration Date: 30 SEP 2016

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: 1201120507104
Free TAG Mean: 15 634
PQ Low Control Mean: 80
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	Quality Control Analyst		16 DEC 2013
Review/Approval	Staci Leffel	Manufacturing Manager		16 DEC 2013
Review/Approval	Richard Mann	Quality Engineer		16 DEC 2013



CERTIFICATE OF ANALYSIS