



CERTIFICATE OF ANALYSIS

COA – 04338

Product Description: MSD® Performance Qualification Kit, Sector Imager

Catalog #/Size: R31QQ-3

Lot Number: PQ300037

Expiration Date: 31 MAR 2016

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

SECTOR Imager Plate Lot Number: 21E0EA
SECTOR Imager Plate Catalog Number: L15XA
SECTOR Imager Plate Expiration Date: 31 MAR 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
PQ Low Control Lot Number: Y0320002
PQ Low Control Catalog Number: R92TE
PQ Low Control Expiration Date: 31 JAN 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: 1200050211221
Free TAG Mean: 16 060
PQ Low Control Mean: 89
Dark Signal Mean: -2

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

	Name	Function	Signature	Date
Preparer	Michelle Gao	Quality Technical Specialist		16APR2014
Review/Approval	Mike Eager	Quality Scientist		16Apr2014