



RFE MEMO

Page 1 of 1

Product Description(s): MSD® SECTOR Imager Performance Qualification Kit

Catalog Number(s) / Size(s): R31QQ-3

Lot Number: PQ300099

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of PQ300099 has been extended to 28 Feb 2019.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
SECTOR Imager Plate	21G0QA	28 FEB 2018	28 FEB 2019

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	21 FEB 2018

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® SECTOR Imager Performance Qualification Kit

Catalog #/Size: R31QQ-3

Lot Number: PQ300099

Expiration Date: 28 FEB 2018

Statement: This product was manufactured and tested according to approved MSD documents.

SECTOR Imager Plate					
Catalog Number	N/A	Lot Number	21G0QA	Expiration Date	28 FEB 2018

Free Tag 15,000					
Catalog Number	N/A	Lot Number	Y0260157	Expiration Date	30 JUN 2020

PQ Low Control					
Catalog Number	R92TA	Lot Number	Y0110066	Expiration Date	29 FEB 2020

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

MSD QC Results:					
Using Instrument	Sector Imager		S/N	1201130115105	
Free TAG Mean	15 573	PQ Low Control Mean	82	Dark Signal Mean	1

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	05 OCT 2017