



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD<sup>®</sup> SECTOR Imager Performance Qualification Kit

**Catalog Number:** R31QQ-3

**Lot Number:** PQ300129

**Expiration Date:** 31 JUL 2021

**Storage Temperature:** Room Temperature

Components	Catalog Number	Lot Number	Expiration Date
SECTOR Imager Plate	L15XA	21J35A	31 JUL 2021
MSD Free Tag ECL 15000	N/A	Y0260163	28 FEB 2022
PQ Low Control	R92TE	Y0110074	31 MAY 2022

Test	QC Obtained Mean Signal	Signal Specifications	
		Minimal Value	Maximal Value
MSD Free Tag ECL 15000	14566	12 000	18 000
PQ Low Control	62	25	100
Dark Noise	1	-16	16
Test	%CV and SD Specification		
MSD Free Tag ECL 15000	0-6.0%		
PQ Low Control SD	0-25		
Dark Noise SD	0-16		

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 OCT 2019