



CERTIFICATE OF ANALYSIS

Product Description: SECTOR Imager Performance Qualification Kit
Catalog Number: R31QQ-3
Lot Number: PQ300124
Expiration Date: 30 Apr 2021

Statement: This product was manufactured and tested according to approved MSD Standard Operating Procedures.

Components for Testing SECTOR Imager Performance Qualification Kit	Catalog Number	Lot Number	Expiration Date
SECTOR Imager Plate	L15XA-3	21J1EA	30 Apr 2021
MSD Free Tag ECL 15000	N/A	Y0260163	28 Feb 2022
PQ Low Control	R92TE-2	Y0110073	31 Jan 2022

SECTOR Imager Performance Qualification Kit Specifications			
Test	QC Obtained Mean Signal	Signal Specifications	
		Minimal Value	Maximal Value
MSD Free Tag ECL 15000	15,922	12,000	18,000
PQ Low Control	80	25	100
Dark Noise	-1	-16	16
Test	%CV and SD Specification		
MSD Free Tag ECL 15000	0 % - 6 %		
PQ Low Control SD	0 - 25		
Dark Noise SD	0 - 16		

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality		16 MAY 2019