



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® SECTOR Imager Performance Qualification Kit

Catalog #/Size: R31QQ-3

Lot Number: PQ300119

Expiration Date: 31 OCT 2020

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

| SECTOR Imager Plate | | | | | |
|---------------------|---------------|------------|--------|-----------------|-------------|
| Catalog Number | L15XA / L11XA | Lot Number | 21I3OA | Expiration Date | 31 OCT 2020 |

| Free Tag 15,000 | | | | | |
|-----------------|-----|------------|----------|-----------------|-------------|
| Catalog Number | N/A | Lot Number | Y0260160 | Expiration Date | 30 NOV 2021 |

| PQ Low Control | | | | | |
|----------------|-------|------------|----------|-----------------|-------------|
| Catalog Number | R92TE | Lot Number | Y0320014 | Expiration Date | 31 MAR 2021 |

| 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 | 15 778 | PQ Low Control Mean | 75 | Dark Signal Mean | -0.9 |

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> | 14 FEB 2019 |