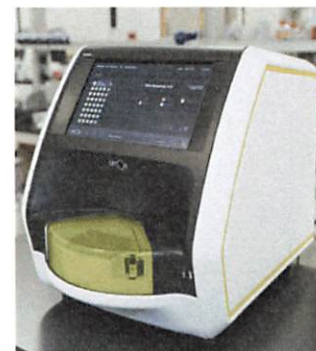


| RQT v2 (5 Plate) | |
|------------------|--------|
| Product Number: | 103642 |
| Lot Number: | N/A |
| Expiration: | N/A |
| Platform(s): | SR-X |

| RQT v2 (1 Plate) | |
|------------------|------------|
| Product Number: | 103641 |
| Lot Number: | 504445 |
| Expiration: | 1-Jan-2025 |
| Platform(s): | SR-X |



Data below represents results generated on the Simoa® SR-X Analyzer.

| Component | Part Number | Lot Number |
|--------------|-------------|------------------|
| RQT v2 Plate | 103891 | 430602 |
| RGP Reagent | 103159 | N/A ¹ |

¹ Reagents are not Kit Lot Specific

| | Acceptance Criteria | Instrument 1 | Instrument 2 |
|----------------------------------|---------------------|--------------|--------------|
| Data Yield | ≥95% | 100 % | 95% |
| Low AEB Avg. | 0.06 - 0.30 | 0.07 | 0.08 |
| High AEB Avg. | 2 - 15 | 9 | 9 |
| Low AEB Number of Beads Avg. | ≥7700 | 22185 | 22689 |
| Low AEB Number of Beads CV | ≤35% CV | 9 % | 9 % |
| High AEB Number of Beads Avg. | ≥7700 | 21629 | 18813 |
| High AEB Number of Beads CV | ≤35% CV | 5 % | 7 % |
| AEB Bead Fill % Avg. | ≥12.5% | 15.9 % | 15.5 % |
| AEB Bead Fill % CV | ≤20% | 6 % | 9 % |
| Isingle | ≥200 | 274 | 341 |
| Multiplex Bead Number (ID: 0001) | ≥1500 beads/plex | 4011 | 2907 |
| Multiplex Bead Number (ID: 0002) | ≥1500 beads/plex | 3298 | 2343 |
| Multiplex Bead Number (ID: 0021) | ≥1500 beads/plex | 3796 | 2633 |
| Multiplex Bead Number (ID: 0201) | ≥1500 beads/plex | 3627 | 2550 |
| Multiplex Bead Number (ID: 2001) | ≥1500 beads/plex | 3856 | 2642 |

Review/Approval

TAM LEE DOC CONTROL II, QA Name Lee 19 NOV 2024
 Name Title Signature/Date