

| 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:      | 504718      |
| Expiration:      | 10-Mar-2026 |
| Platform(s):     | SR-X        |



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

| Component    | Part Number | Lot Number       |
|--------------|-------------|------------------|
| RQT v2 Plate | 103891      | 525301           |
| RGP Reagent  | 103159      | N/A <sup>1</sup> |

<sup>1</sup> Reagents are not Kit Lot Specific

|                                  | Acceptance Criteria | Instrument 1 | Instrument 2 |
|----------------------------------|---------------------|--------------|--------------|
| Data Yield                       | ≥95%                | 99 %         | 99 %         |
| Low AEB Avg.                     | 0.06 - 0.30         | 0.17         | 0.17         |
| High AEB Avg.                    | 2 - 15              | 7            | 6            |
| Low AEB Number of Beads Avg.     | ≥7700               | 22207        | 23113        |
| Low AEB Number of Beads CV       | ≤35% CV             | 8 %          | 9 %          |
| High AEB Number of Beads Avg.    | ≥7700               | 22185        | 24012        |
| High AEB Number of Beads CV      | ≤35% CV             | 8 %          | 11 %         |
| AEB Bead Fill % Avg.             | ≥12.5%              | 16.2 %       | 16.8 %       |
| AEB Bead Fill % CV               | ≤20%                | 7 %          | 9 %          |
| Isingle                          | ≥200                | 287          | 247          |
| Multiplex Bead Number (ID: 0001) | ≥1500 beads/plex    | 2712         | 2855         |
| Multiplex Bead Number (ID: 0002) | ≥1500 beads/plex    | 3297         | 3422         |
| Multiplex Bead Number (ID: 0021) | ≥1500 beads/plex    | 3740         | 3888         |
| Multiplex Bead Number (ID: 0201) | ≥1500 beads/plex    | 3680         | 3766         |
| Multiplex Bead Number (ID: 2001) | ≥1500 beads/plex    | 3676         | 3750         |

Review/Approval

*Dhruvi Patel*

Name

*QA Specialist II, QA*

Title

*DP/18 Sep 2025*

Signature/Date