

RQT v2 (5 Plate)	
Product Number:	103642
Lot Number:	504613
Expiration:	30-Oct-2025
Platform(s):	SR-X

RQT v2 (1 Plate)	
Product Number:	103641
Lot Number:	504591
Expiration:	30-Oct-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	512001
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	100 %	98 %
Low AEB Avg.	0.06 - 0.30	0.20	0.19
High AEB Avg.	2 - 15	4	4
Low AEB Number of Beads Avg.	≥7700	20886	21970
Low AEB Number of Beads CV	≤35% CV	18 %	6 %
High AEB Number of Beads Avg.	≥7700	20936	22605
High AEB Number of Beads CV	≤35% CV	10 %	8 %
AEB Bead Fill % Avg.	≥12.5%	15.9 %	15.7 %
AEB Bead Fill % CV	≤20%	14 %	7 %
Isingle	≥200	297	210
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	4306	3385
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	3448	3195
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3890	3631
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3354	3032
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	3454	3224

Review/Approval

Dhruvi Patel

Name

Doc Control II, QA

Title

DP 28 May 2025

Signature/Date