

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:	536622
Expiration:	6-May-2026
Platform(s):	SR-X

Component	Part Number	Lot Number
RQT v2 Plate	103891	531001
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

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

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	97%	98%
Low AEB Avg.	0.06 - 0.30	0.16	0.14
High AEB Avg.	2 - 15	8	8
Low AEB Number of Beads Avg.	≥7700	23534	23680
Low AEB Number of Beads CV	≤35% CV	8%	12%
High AEB Number of Beads Avg.	≥7700	23569	24443
High AEB Number of Beads CV	≤35% CV	9%	16%
AEB Bead Fill % Avg.	≥12.5%	17.3%	17.4%
AEB Bead Fill % CV	≤20%	7%	11%
Isingle	≥200	265	228
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	3123	3236
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	3478	3688
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3439	3631
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	4098	4279
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	3979	4157

Review/Approval

Dhvani Patel QA Specialist II DP/20 Nov 2025
Name Title Signature/Date