

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:	504456
Expiration:	21-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	432603
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	98%	99%
Low AEB Avg.	0.06 - 0.30	0.10	0.11
High AEB Avg.	2 - 15	9	9
Low AEB Number of Beads Avg.	≥7700	21631	23322
Low AEB Number of Beads CV	≤35% CV	13 %	7 %
High AEB Number of Beads Avg.	≥7700	21815	20833
High AEB Number of Beads CV	≤35% CV	8 %	6 %
AEB Bead Fill % Avg.	≥12.5%	15.4 %	16.4 %
AEB Bead Fill % CV	≤20%	10 %	6 %
Isingle	≥200	240	300
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	3164	2629
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	2812	2515
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3250	2861
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3115	2796
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	3271	2869

Review/Approval

Dipesh Patel
Name

Documentation Lead
Title

06 DEC 2024
Signature/Date