

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:	504409
Expiration:	16-Dec-2024
Platform(s):	SR-X



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

Component	Part Number	Lot Number
RQT v2 Plate	103891	429004
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%	97%
Low AEB Avg.	0.06 - 0.30	0.14	0.17
High AEB Avg.	2 - 15	9	10
Low AEB Number of Beads Avg.	≥7700	23101	23634
Low AEB Number of Beads CV	≤35% CV	9%	9%
High AEB Number of Beads Avg.	≥7700	21642	20061
High AEB Number of Beads CV	≤35% CV	7%	5%
AEB Bead Fill % Avg.	≥12.5%	16.1%	16.5%
AEB Bead Fill % CV	≤20%	8%	7%
Isingle	≥200	286	328
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	4123	3733
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	3158	3258
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3583	3722
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3742	3918
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	3589	3707

Review/Approval

*Dhruvi Patel*    *Doc Control I, QA*    *DP 23 Oct 2024*

Name

Title

Signature/Date