

RQT v2 (5 Plate)	
Product Number:	103642
Lot Number:	504096
Expiration:	22-Jun-2024
Platform(s):	SR-X

RQT v2 (1 Plate)	
Product Number:	103641
Lot Number:	N/A
Expiration:	N/A
Platform(s):	SR-X



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

Component	Part Number	Lot Number
RQT v2 Plate	103891	335601
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%	96%
Bead Fill % Avg.	>12.5%	15.8%	12.7%
Bead Fill % CV	<20%	8%	11%
Isingle	>200	292	264
Low AEB Range	0.06 - 0.30	0.15	0.15
High AEB Range	2 - 15	9	10
High Bead Number	>7700	22926	17637
High Bead CV	<35% CV	8%	16%
Low AEB Bead Number	>7700	21346	15547
Low AEB Bead CV	<35% CV	7%	13%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	4048	3882
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	3328	2266
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	3440	2359
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	3867	2639
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	3603	2448

Review/Approval

Cody Carlson QC Manager  
 Name Title Signature/Date  
 [Signature] 27 Dec 2023