

RQT v2 (5 Plate)	
Product Number:	103642
Lot Number:	504126
Expiration:	1-Aug-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	403204
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	>95%	98%	99%
Bead Fill % Avg.	>12.5%	17.4%	17.6%
Bead Fill % CV	<20%	6%	7%
Isingle	>200	287	328
Low AEB Range	0.06 - 0.30	0.14	0.16
High AEB Range	2 - 15	8	9
High Bead Number	>7700	22907	24525
High Bead CV	<35% CV	7%	8%
Low AEB Bead Number	>7700	23597	24753
Low AEB Bead CV	<35% CV	7%	6%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	3655	3548
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	3369	3410
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	3649	3705
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	3886	4128
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	3562	3670

Review/Approval

Cody Carlson
Name
QC Manager
Title
Cody Carlson 05Feb2024
Signature/Date