

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:	504343
Expiration:	12-Jun-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	410307
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	99 %	98 %
Low AEB Avg.	0.06 - 0.30	0.14	0.12
High AEB Avg.	2 - 15	7	6
Low AEB Number of Beads Avg.	≥7700	19389	18561
Low AEB Number of Beads CV	≤35% CV	11 %	13 %
High AEB Number of Beads Avg.	≥7700	18600	19360
High AEB Number of Beads CV	≤35% CV	17 %	12 %
AEB Bead Fill % Avg.	≥12.5%	14.5 %	14.7 %
AEB Bead Fill % CV	≤20%	12 %	8 %
Isingle	≥200	304	265
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	4465	3970
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	2650	3092
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	2667	3084
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3051	3459
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	2939	3371

Review/Approval

Cody Carlson
Name

QC Manager
Title

[Signature] 07 AUG 2024
Signature/Date