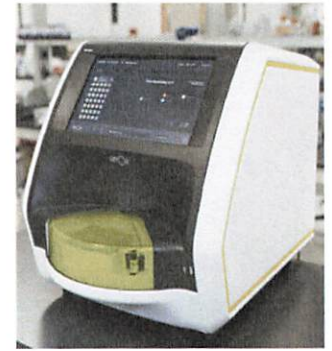


RQT v2 (5 Plate)	
Product Number:	103642
Lot Number:	504237
Expiration:	29-Jul-2024
Platform(s):	SR-X

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	415204
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	0.13	0.10
High AEB Avg.	2 - 15	8	8
Low AEB Number of Beads Avg.	≥7700	19985	20666
Low AEB Number of Beads CV	≤35% CV	12 %	8 %
High AEB Number of Beads Avg.	≥7700	19719	21042
High AEB Number of Beads CV	≤35% CV	6 %	9 %
AEB Bead Fill % Avg.	≥12.5%	14.7 %	15.3 %
AEB Bead Fill % CV	≤20%	9 %	7 %
Isingle	≥200	328	253
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	3028	3778
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	2711	2412
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	2903	2589
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3330	3101
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	2923	2645

Review/Approval

Cody Carlson

QC Manager

[Signature] 06 JUN 2024

Name

Title

Signature/Date