

RQT v2 (5 Plate)	
Product Number:	103642
Lot Number:	504187
Expiration:	29-Aug-2025
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	408201
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	98 %	98 %
Low AEB Avg.	0.06 - 0.30	0.15	0.14
High AEB Avg.	2 - 15	6	5
Low AEB Number of Beads Avg.	≥7700	31442	19918
Low AEB Number of Beads CV	≤35% CV	13%	14 %
High AEB Number of Beads Avg.	≥7700	28646	20161
High AEB Number of Beads CV	≤35% CV	22 %	10 %
AEB Bead Fill % Avg.	≥12.5%	21.6%	15.2 %
AEB Bead Fill % CV	≤20%	16 %	11 %
Isingle	≥200	381	292
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	3903	3231
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	3808	2643
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3758	2710
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3934	3007
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	4317	2719

Review/Approval

Cody Carlson

QC Manager

Cody Carlson 25 MAR 2024
Signature/Date

Name

Title