

RQT v2 (5 Plate)	
Product Number:	103642
Lot Number:	504185
Expiration:	15-Sep-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	407501
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥ 95%	99 %	98 %
Bead Fill % Avg.	≥ 12.5%	19.3 %	16.3 %
Bead Fill % CV	≤ 20%	19 %	8 %
Isingle	≥ 200	340	248
Low AEB Avg.	0.06 - 0.30	0.15	0.14
High AEB Avg.	2 - 15	7	8
High Bead Number	≥ 7700	23289	21398
High AEB Bead CV	≤ 35% CV	21 %	9 %
Low AEB Bead Number	≥ 7700	27674	21719
Low AEB Bead CV	≤ 35% CV	19 %	10 %
Multiplex Bead Number (ID: 0001)	≥ 1500 beads/plex	4743	4008
Multiplex Bead Number (ID: 0002)	≥ 1500 beads/plex	3848	3108
Multiplex Bead Number (ID: 0021)	≥ 1500 beads/plex	4070	3254
Multiplex Bead Number (ID: 0201)	≥ 1500 beads/plex	4610	3656
Multiplex Bead Number (ID: 2001)	≥ 1500 beads/plex	4416	3563

Review/Approval

Cody Carlson

QC Manager

Cody Carlson 20 MAR 2024

Name

Title

Signature/Date