

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:	504112
Expiration:	14-Aug-2024
Platform(s):	SR-X

Component	Part Number	Lot Number
RQT v2 Plate	103891	404503
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific



Data below represents results generated on the Simoa® SR-X Analyzer.

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	>95%	97%	99%
Bead Fill % Avg.	>12.5%	16.8%	14.3%
Bead Fill % CV	<20%	17%	13%
Isingle	>200	377	278
Low AEB Range	0.06 - 0.30	0.14	0.14
High AEB Range	2 - 15	7	7
High Bead Number	>7700	20620	16506
High Bead CV	<35% CV	20%	13%
Low AEB Bead Number	>7700	25054	20946
Low AEB Bead CV	<35% CV	13%	7%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	4345	3959
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	3523	2966
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	3683	3139
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	4053	3560
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	3758	3211

Review/Approval

Cody Carlson
Name

QC Manager
Title

[Signature] 20 FEB 2024
Signature/Date