

RQT v2 (5 Plate)	
Product Number:	103642
Lot Number:	504082
Expiration:	14-Jun-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	334802
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	>95%	100%	100%
Bead Fill % Avg.	>12.5%	16.2%	15.2%
Bead Fill % CV	<20%	10%	7%
Isingle	>200	357	335
Low AEB Range	0.06 - 0.30	0.17	0.17
High AEB Range	2 - 15	11	11
High Bead Number	>7700	22127	20415
High Bead CV	<35% CV	9%	7%
Low AEB Bead Number	>7700	22565	21265
Low AEB Bead CV	<35% CV	11%	8%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	4288	5114
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	3544	3393
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	3438	3276
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	4235	4077
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	3712	3551

Review/Approval

ROBERT SOMOGYI

QC MANAGER

Signature/Date

Name

Title

Signature/Date