

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:	504125
Expiration:	1-Aug-2024
Platform(s):	SR-X



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

Component	Part Number	Lot Number
RQT v2 Plate	103891	403203
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	>95%	96%	96%
Bead Fill % Avg.	>12.5%	21.5%	15.7%
Bead Fill % CV	<20%	13%	10%
Isingle	>200	317	248
Low AEB Range	0.06 - 0.30	0.16	0.13
High AEB Range	2 - 15	9	8
High Bead Number	>7700	28531	20430
High Bead CV	<35% CV	16%	16%
Low AEB Bead Number	>7700	29393	19865
Low AEB Bead CV	<35% CV	12%	13%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	4229	4278
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	3943	2524
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	4322	2742
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	4959	3118
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	4249	2731

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

[Signature] 02 FEB 2024
Signature/Date