

Product Number:	103642
Lot Number:	504029
Expiration:	25-Apr-2024
Platform(s):	Simoa SR-X Analyzer



Component	Item Number	Lot Number
RQT v2 Plate	103891	329801
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

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

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	>95%	98 %	99 %
Bead Fill % Avg.	>12.5%	13.5 %	15.1 %
Bead Fill % CV	<20%	18 %	14 %
Isingle	>200	274	322
Low AEB Range	0.06 - 0.30	0.15	0.15
High AEB Range	2 - 15	7	7
High Bead Number	>7700	14504	15847
High Bead CV	<35% CV	15 %	15 %
Low AEB Bead Number	>7700	21275	22543
Low AEB Bead CV	<35% CV	6 %	9 %
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	4163	4570
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	2432	3274
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	2518	3321
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	2841	3721
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	2513	3300

Review/Approval

ROBERT SMOGYI
Name

QC MANAGER
Title

[Signature] 31 OCT 2023
Signature/Date