

Product Number:	103642
Lot Number:	503971
Expiration:	11-Mar-2024
Platform(s):	Simoa SR-X Analyzer



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

Component	Item Number	Lot Number
RQT v2 Plate	103891	325405
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	>95%	98 %	100 %
Bead Fill % Avg.	>12.5%	15.3 %	15.7 %
Bead Fill % CV	<20%	7 %	9 %
Isingle	>200	374	253
Low AEB Range	0.06 - 0.30	0.15	0.12
High AEB Range	2 - 15	6	6
High Bead Number	>7700	19731	20422
High Bead CV	<35% CV	10 %	14 %
Low AEB Bead Number	>7700	21106	21664
Low AEB Bead CV	<35% CV	8 %	8 %
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	4033	4813
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	2591	2752
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	2883	3013
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	3046	3139
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	2721	2852

Review/Approval

ROBERT SOMOGYI QC, MANAGER
Name Title

12 SEP 2023
Signature/Date