

Product Number:	103642
Lot Number:	504048
Expiration:	10-May-2024
Platform(s):	Simoa SR-X Analyzer



Component	Item Number	Lot Number
RQT v2 Plate	103891	331402
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%	96%	100%
Bead Fill % Avg.	>12.5%	20.4%	15.2%
Bead Fill % CV	<20%	18%	11%
Isingle	>200	385	283
Low AEB Range	0.06 - 0.30	0.17	0.16
High AEB Range	2 - 15	8	8
High Bead Number	>7700	27057	20645
High Bead CV	<35% CV	25%	11%
Low AEB Bead Number	>7700	28746	19978
Low AEB Bead CV	<35% CV	13%	12%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	4647	4566
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	4119	2819
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	4577	3120
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	5117	3444
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	4617	3164

Review/Approval

Sania Khillan
Name

Senior QC Analyst
Title

Sania Khillan 14/NOV/2023
Signature/Date