

Product Number:	103642
Lot Number:	503830
Expiration:	16-Sep-2023
Platform(s):	Simoa SR-X Analyzer



Component	Item Number	Lot Number
RQT v2 Plate	103891	307501
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%	99%	98%
Bead Fill % Avg.	>12.5%	14.5%	14.5%
Bead Fill % CV	<20%	9%	9%
Isingle	>250	263	293
Low AEB Range	0.06 - 0.30	0.13	0.15
High AEB Range	2 - 15	10	11
High Bead Number	>7700	18710	18638
High Bead CV	<35% CV	14%	12%
Low AEB Bead Number	>7700	20851	20387
Low AEB Bead CV	<35% CV	8%	10%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	4015	3938
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	2692	2854
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	2788	2977
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	3325	3452
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	2972	3127

Review/Approval

Sonia Khillan
Name

QC

Title

Sonia Khillan 07 APR 2023
Signature/Date