

Product Number:	103642
Lot Number:	503888
Expiration:	24-Nov-2023
Platform(s):	Simoa SR-X Analyzer



Component	Item Number	Lot Number
RQT v2 Plate	103891	314401 ¹
RGP Reagent	103159	N/A ²

¹ RQT v2 Plate control Lot# 314401A

² 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%	97%	99%
Bead Fill % Avg.	>12.5%	21.6%	18.6%
Bead Fill % CV	<20%	14%	12%
Isingle	>200	288	242
Low AEB Range	0.06 - 0.30	0.13	0.11
High AEB Range	2 - 15	11	9
High Bead Number	>7700	25952	123888
High Bead CV	<35% CV	19%	16%
Low AEB Bead Number	>7700	28865	26140
Low AEB Bead CV	<35% CV	16%	10%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	5392	5879
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	3832	3418
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	3949	3464
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	4733	4127
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	4222	3718

Review/Approval

Sonia Khillan
Name

QC

Title

Sonia 12/1/23 06 JUN 2023
Signature/Date