

Product Number:	103642
Lot Number:	503937
Expiration:	7-Feb-2024
Platform(s):	Simoa SR-X Analyzer



Component	Item Number	Lot Number
RQT v2 Plate	103891	321904
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%	98%
Bead Fill % Avg.	>12.5%	15.5%	13.7%
Bead Fill % CV	<20%	16%	17%
Isingle	>200	255	289
Low AEB Range	0.06 - 0.30	0.12	0.14
High AEB Range	2 - 15	6	6
High Bead Number	>7700	19365	13894
High Bead CV	<35% CV	13%	28%
Low AEB Bead Number	>7700	21544	19542
Low AEB Bead CV	<35% CV	23%	13%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	3711	3795
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	2321	1856
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	2365	1871
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	2765	2245
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	2671	2162

Review/Approval

ROBERT SANDOZY QC, MANAGER
Name Title

[Signature] 15 AUG 2023
Signature/Date

Quanterix Corporation

900 Middlesex Turnpike, Building 1, Billerica, MA 01821

techsupport@quanterix.com

©Quanterix Corp. Simoa® is a registered trademark of Quanterix Corp.

For research use only. Not for use in diagnostic procedures