

Product Number:	103642
Lot Number:	503441
Expiration:	27 Oct 2022
Platform(s):	Simoa SR-X Analyzer



Component	Item Number	Lot Number
RQT v2 Plate	103891	211705
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%	100%
Bead Fill % Avg.	>12.5%	15.4%	15.0%
Bead Fill % CV	<20%	5.55%	10.4%
Isingle	>250	344	282
Low AEB Range	0.06 - 0.30	0.100	0.100
High AEB Range	2 - 15	8.78	8.50
Low AEB Bead Number	>7700	21581	19317
High AEB Bead Number	>7700	21028	18068
Low AEB Bead CV	<35% CV	6.6%	14.3%
High AEB Bead CV	<35% CV	6.2%	23.4%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	5029	2975
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	3031	3124
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	2710	2484
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	3222	2530
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	3049	2793

Review/Approval

Matt Smart

Director, Quality Control

MA 23 May 2022

Name

Title

Signature/Date

Quanterix Corporation

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

900 Middlesex Turnpike, Building 1, Billerica, MA 01821

For research use only. Not for use in diagnostic procedures

techsupport@quantexix.com