

Product Number:	103642
Lot Number:	503809
Expiration:	10-Aug-2023
Platform(s):	Simoa SR-X Analyzer



Component	Item Number	Lot Number
RQT v2 Plate	103891	304105
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%	17.5 %	17.1 %
Bead Fill % CV	<20%	10 %	9.0 %
Isingle	>250	292	291
Low AEB Range	0.06 - 0.30	0.15	0.16
High AEB Range	2 - 15	12.8	12.8
High and Low AEB Bead Number	>7700	23408	22185
High and Low AEB Bead CV	<35% CV	9.5 %	7.5 %
Multiplex Bead Number	>1500 beads/plex	3738	3392

Review/Approval

Sonia Khiln
Name

QC

Title

Sonia Khiln 1/15 FEB 2023
Signature/Date