

Product Number: 103642
Lot Number: 503908
Expiration: 23-Dec-2023
Platform(s): Simoa SR-X Analyzer



Data below
represents
results
generated on
the Simoa® SR-
X Analyzer.

Component	Item Number	Lot Number	Control Lot Number
RQT v2 Plate	103891	317402	317402A
RGP Reagent	103159	N/A ¹	N/A ¹

¹ RGP is not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	>95%	98%	98%
Bead Fill % Avg.	>12.5%	14.9%	13.4%
Bead Fill % CV	<20%	12%	17%
Isingle	>250	381	340
Low AEB Range	0.06 - 0.30	0.16	0.16
High AEB Range	2 - 15	6.82	6.30
High Bead Number	>7700	17319	14682
High Bead CV	<35% CV	12%	18%
Low AEB Bead Number	>7700	20793	19635
Low AEB Bead CV	<35% CV	12%	17%
Multiplex Bead Number (ID: 0001)	>1500 beads/plex	3630	5655
Multiplex Bead Number (ID: 0002)	>1500 beads/plex	1888	1931
Multiplex Bead Number (ID: 0021)	>1500 beads/plex	2234	2248
Multiplex Bead Number (ID: 0201)	>1500 beads/plex	2652	2676
Multiplex Bead Number (ID: 2001)	>1500 beads/plex	2217	2257

Review/Approval

Sonia Khilkin
Name

QC Manager
Title

Soni Koll 13 Jul 2023
Signature/Date