

RQT v2 (5 Plate)	
Product Number:	103642
Lot Number:	N/A
Expiration:	N/A
Platform(s):	SR-X

RQT v2 (1 Plate)	
Product Number:	103641
Lot Number:	504445
Expiration:	1-Feb-2025
Platform(s):	SR-X



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

Component	Part Number	Lot Number
RQT v2 Plate	103891	430602
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	100 %	95%
Low AEB Avg.	0.06 - 0.30	0.07	0.08
High AEB Avg.	2 - 15	9	9
Low AEB Number of Beads Avg.	≥7700	22185	22689
Low AEB Number of Beads CV	≤35% CV	9 %	9 %
High AEB Number of Beads Avg.	≥7700	21629	18813
High AEB Number of Beads CV	≤35% CV	5 %	7 %
AEB Bead Fill % Avg.	≥12.5%	15.9 %	15.5 %
AEB Bead Fill % CV	≤20%	6 %	9 %
Isingle	≥200	274	341
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	4011	2907
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	3298	2343
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3796	2633
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3627	2550
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	3856	2642

Review/Approval

Dipesh Patel
Name

Doc Control II QA
Title

23 JAN 2025
Signature/Date