

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:	504545
Expiration:	19-Jun-2025
Platform(s):	SR-X



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

Component	Part Number	Lot Number	Control Lot Number
RQT v2 Plate	103891	507802	507802A
RGP Reagent	103159	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	100 %	96 %
Low AEB Avg.	0.06 - 0.30	0.20	0.23
High AEB Avg.	2 - 15	4	5
Low AEB Number of Beads Avg.	≥7700	21987	23756
Low AEB Number of Beads CV	≤35% CV	19 %	15 %
High AEB Number of Beads Avg.	≥7700	22476	23340
High AEB Number of Beads CV	≤35% CV	7 %	19 %
AEB Bead Fill % Avg.	≥12.5%	16.0 %	17.0 %
AEB Bead Fill % CV	≤20%	14 %	16 %
Isingle	≥200	216	284
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	3186	4048
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	3117	2643
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3406	2976
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3719	3275
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	3709	3277

Review/Approval

Dhruv Patel

DOC Control, II QA

DP 15 Apr 2025

Name

Title

Signature/Date