

RQT v2 (5 Plate)	
Product Number:	103642
Lot Number:	504697
Expiration:	14-Feb-2026
Platform(s):	SR-X

RQT v2 (1 Plate)	
Product Number:	103641
Lot Number:	504693
Expiration:	14-Feb-2026
Platform(s):	SR-X



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

Component	Part Number	Lot Number
RQT v2 Plate	103891	522604
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	100 %	99 %
Low AEB Avg.	0.06 - 0.30	0.14	0.14
High AEB Avg.	2 - 15	8	8
Low AEB Number of Beads Avg.	≥7700	23339	21548
Low AEB Number of Beads CV	≤35% CV	11 %	15 %
High AEB Number of Beads Avg.	≥7700	22237	21379
High AEB Number of Beads CV	≤35% CV	21 %	16 %
AEB Bead Fill % Avg.	≥12.5%	16.5%	15.6 %
AEB Bead Fill % CV	≤20%	16 %	14 %
Isingle	≥200	267	225
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	3215	2927
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	3278	3093
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3633	3467
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3903	3700
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	3474	3263

Review/Approval

Dhruvi Patel

QA Specialist II, QA

SDR | *21 Aug 2025*

Name

Title

Signature/Date