

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:	504529
Expiration:	13-Jun-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	507208
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	100 %	98 %
Low AEB Avg.	0.06 - 0.30	0.09	0.08
High AEB Avg.	2 - 15	5	5
Low AEB Number of Beads Avg.	≥7700	21705	20957
Low AEB Number of Beads CV	≤35% CV	9 %	12 %
High AEB Number of Beads Avg.	≥7700	22991	23097
High AEB Number of Beads CV	≤35% CV	8 %	8 %
AEB Bead Fill % Avg.	≥12.5%	16.3%	16.2 %
AEB Bead Fill % CV	≤20%	4 %	7 %
Isingle	≥200	286	245
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	3268	2991
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	3563	3324
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3635	3359
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	4046	3758
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	3781	3532

Review/Approval

Dipesh Patel
Name

Str. Doc Control, QA
Title

20 MAR 2025
Signature/Date