

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:	504361
Expiration:	23-Oct-2024
Platform(s):	SR-X



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

Component	Part Number	Lot Number
RQT v2 Plate	103891	423606
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	98 %	96 %
Low AEB Avg.	0.06 - 0.30	0.06	0.07
High AEB Avg.	2 - 15	4	5
Low AEB Number of Beads Avg.	≥7700	19858	22427
Low AEB Number of Beads CV	≤35% CV	16 %	7 %
High AEB Number of Beads Avg.	≥7700	20393	21825
High AEB Number of Beads CV	≤35% CV	15 %	15 %
AEB Bead Fill % Avg.	≥12.5%	14.5%	16.0 %
AEB Bead Fill % CV	≤20%	15 %	11 %
Isingle	≥200	234	291
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	3199	3808
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	2881	3488
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3072	3670
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3144	3770
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	3212	3810

Review/Approval

Dhruvi Patel *Doc Control I, QA* *DP 20 Sep 2024*
 Name Title Signature/Date