

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:	504351
Expiration:	15-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	422803
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

	Acceptance Criteria	Instrument 1	Instrument 2
Data Yield	≥95%	100%	100%
Low AEB Avg.	0.06 - 0.30	0.07	0.07
High AEB Avg.	2 - 15	5	4
Low AEB Number of Beads Avg.	≥7700	22224	23012
Low AEB Number of Beads CV	≤35% CV	9%	10%
High AEB Number of Beads Avg.	≥7700	17940	18282
High AEB Number of Beads CV	≤35% CV	9%	14%
AEB Bead Fill % Avg.	≥12.5%	15.2%	16.4%
AEB Bead Fill % CV	≤20%	11%	11%
Isingle	≥200	263	227
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	2821	3328
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	2786	3119
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	3202	3431
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	3277	3742
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	3216	3583

Review/Approval

Difesh Patel
Name

Doc Control - 2 QA
Title

mt / 26 Aug 2024
Signature/Date