

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:	504492
Expiration:	23-Apr-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	502302	502302A
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%	95%	99 %
Low AEB Avg.	0.06 - 0.30	0.06	0.07
High AEB Avg.	2 - 15	8	9
Low AEB Number of Beads Avg.	≥7700	32288	21085
Low AEB Number of Beads CV	≤35% CV	11%	12%
High AEB Number of Beads Avg.	≥7700	30462	19056
High AEB Number of Beads CV	≤35% CV	10%	12%
AEB Bead Fill % Avg.	≥12.5%	22.7%	14.9%
AEB Bead Fill % CV	≤20%	10%	10%
Isingle	≥200	230	248
Multiplex Bead Number (ID: 0001)	≥1500 beads/plex	4966	3804
Multiplex Bead Number (ID: 0002)	≥1500 beads/plex	4506	3317
Multiplex Bead Number (ID: 0021)	≥1500 beads/plex	4943	3307
Multiplex Bead Number (ID: 0201)	≥1500 beads/plex	5027	3681
Multiplex Bead Number (ID: 2001)	≥1500 beads/plex	5206	3570

Review/Approval

Direk Patel Doc control / IT QA et 03 Feb 2025
Name Title Signature/Date