

Product Number:	103187
Lot Number:	503219
Expiration:	28-Feb-2022

Component	Item Number	Lot Number
RQT Plate	103095	124201
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Means of Analysis	Result
Simoa SR-X Analyzer	PASS

Review/Approval Shivani Goel QC Supervisor Shivani Goel 16 Nov 2021
 Name Title Signature/Date