

Product Number:	103187
Lot Number:	502806
Expiration:	14-Sep-2021

Component	Item Number	Lot Number
RQT Plate	103095	107305
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Means of Analysis	Result
Simoa SR-X Analyzer	PASS

Review/Approval

Junpil Park

QC Supervisor

 19 Mar 2021

Name

Title

Signature/Date