

Product Number:	103187
Lot Number:	502979
Expiration:	7-Dec-2021

Component	Item Number	Lot Number
RQT Plate	103095	115813
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

 17 Jun 21

Name

Title

Signature/Date