

Product Number:	103187
Lot Number:	502727
Expiration:	4-Aug-2021

Component	Item Number	Lot Number
RQT Plate	103095	103501
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
Name

QC Supervisor
Title

 10 Feb 2021
Signature/Date