

Product Number:	103187
Lot Number:	502653
Expiration:	8-Jun-2021

Component	Item Number	Lot Number
RQT Plate	103095	034304
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

 16 Dec 2020
Signature/Date