

Product Number:	103187
Lot Number:	503325
Expiration:	16-Aug-2022

Component	Item Number	Lot Number
RQT Plate	103095	204706
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 Manager

Shivani Goel 22Feb2022

Name

Title

Signature/Date