

Product Number:	103187
Lot Number:	503278
Expiration:	18-Jun-2022

Component	Item Number	Lot Number
RQT Plate	103095	135207
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
Name

QC

Title

Shivani Goel 29 Dec 2021
Signature/Date