

Product Number:	103187
Lot Number:	502543
Expiration:	25-Feb-2021

Component	Item Number	Lot Number
RQT Plate	103095	023804
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 24 Sep 2020

Name

Title

Signature/Date