

Product Number:	103187
Lot Number:	502155
Expiration:	6-Jun-2020

Component	Item Number	Lot Number
RQT Plate	103095	934020
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Means of Analysis	Result
Simoa SR-X Analyzer	PASS

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

12 Dec 2019

Name

Title

Signature/Date