

Product Number:	103187
Lot Number:	502278
Expiration:	6-Sep-2020

Component	Item Number	Lot Number
RQT Plate	103095	006608
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
Name

Assoc. Director, QC
Title

P. Stadler

Signature/Date

10 Mar 2020