

Product Number:	103187
Lot Number:	502235
Expiration:	7-Aug-2020

Component	Item Number	Lot Number
RQT Plate	103095	003801
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 Feb 2020

Name

Title

Signature/Date