

Product Number:	103187
Lot Number:	502060
Expiration:	3-Apr-2020

Component	Item Number	Lot Number
RQT Plate	103095	927604
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Means of Analysis	Result
Simoa SR-X Analyzer	PASS

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*P. Stadler*

15 Oct 2019

Name

Title

Signature/Date