

Product Number:	103187
Lot Number:	501923
Expiration:	1-Jan-20

Component	Item Number	Lot Number
RQT Plate	103095	501910
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*

12 Aug 2019

Name

Title

Signature/Date