

<b>Product Number:</b>	103187
<b>Lot Number:</b>	502456
<b>Expiration:</b>	29-Dec-2020

Component	Item Number	Lot Number
RQT Plate	103095	018103
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

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 15 Jul 2020  
Signature/Date