

<b>Product Number:</b>	103187
<b>Lot Number:</b>	502937
<b>Expiration:</b>	12-Nov-2021

Component	Item Number	Lot Number
RQT Plate	103095	113201
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

Junpil Park  
Name

QC Supervisor  
Title

 27 May 2021  
Signature/Date