

Product Number:	103187
Lot Number:	502696
Expiration:	20-Jul-2021

Component	Item Number	Lot Number
RQT Plate	103095	102010
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

QC Supervisor



21 Jan 2021

Name

Title

Signature/Date