

Product Number:	103187
Lot Number:	503654
Expiration:	01-Feb-2023

Component	Item Number	Lot Number
RQT Plate	103095	221301
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Means of Analysis	Result
Simoa SR-X Analyzer	PASS



Data below represents results generated on the Simoa® SR-X Analyzer.

Review/Approval	<u> Matt Koelaris </u>	<u> QC Supervisor </u>	<u> [Signature] </u>	<u> 23 Aug 2022 </u>
	Name	Title	Signature/Date	