

Product Number:	103187
Lot Number:	503708
Expiration:	17-Apr-2023

Component	Item Number	Lot Number
RQT Plate	103095	229001
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

Matt Kalouris

Name

QC supervisor

Title

[Signature]

Signature/Date

31 Oct 2022