

Product Number:	103187
Lot Number:	503440
Expiration:	1-Nov-2022

Component	Item Number	Lot Number
RQT Plate	103095	212101
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

Matthew Smart

QC Director

[Handwritten Signature] 27 May 2022

Name

Title

Signature/Date