

Product Number:	103187
Lot Number:	502012
Expiration:	4-Mar-2020

Component	Item Number	Lot Number
RQT Plate	103095	924706
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Means of Analysis	Result
Simoa SR-X Analyzer	PASS

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

p. Stadler
Signature/Date

06 Sep 2019