

|                 |             |
|-----------------|-------------|
| Product Number: | 103187      |
| Lot Number:     | 503485      |
| Expiration:     | 23-Dec-2022 |

| Component   | Item Number | Lot Number       |
|-------------|-------------|------------------|
| RQT Plate   | 103095      | 217405           |
| 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   |



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

Review/Approval

Matt Smart      Director, QC      MAT 28 Jan 2022  
 Name                      Title                      Signature/Date