

|                 |             |
|-----------------|-------------|
| Product Number: | 101736      |
| Lot Number:     | 502487      |
| Expiration:     | 27-May-2021 |
| Storage:        | 2-8°C       |

| Component   | Part Number | Lot Number |
|-------------|-------------|------------|
| RGP Reagent | 103159      | 01480      |

| Means of Analysis | PASS/FAIL |
|-------------------|-----------|
| HD-1/HD-X         | Pass      |

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 03 Aug 2020  
Signature/Date