

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 101736      |
| <b>Lot Number:</b>     | 502662      |
| <b>Expiration:</b>     | 20-Nov-2021 |
| <b>Storage:</b>        | 2-8°C       |

| <b>Component</b> | <b>Part Number</b> | <b>Lot Number</b> |
|------------------|--------------------|-------------------|
| RGP Reagent      | 103159             | 03250             |

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

Review/Approval

Junpil Park  
Name

QC, Supervisor  
Title

 29 Dec 2020  
Signature/Date