

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 103591      |
| <b>Lot Number:</b>     | 503413      |
| <b>Expiration:</b>     | 23-Mar-2023 |

| Component        | Part Number | Lot Number       |
|------------------|-------------|------------------|
| Bead Reagent     | 103581      | 209008           |
| Detector Reagent | 103583      | 208808           |
| SBG Reagent      | 103585      | 208703           |
| Control 1        | 103589      | 209101           |
| Control 2        | 103590      | 209101           |
| RGP Reagent      | 103159      | N/A <sup>1</sup> |

<sup>1</sup> RGP is not Kit Lot Specific

|                       |                        |
|-----------------------|------------------------|
| Bead Count (beads/mL) | 1.99 x 10 <sup>7</sup> |
| Bead % Monomeric      | 93%                    |
| Isingle               | 147                    |

**Review/Approval**

Richa Shrestha

QC Supervisor

**Name**

**Title**

**Signature/Date**

*Richa Shrestha* 26APR2022