

<b>Product Number:</b>	101537
<b>Lot Number:</b>	502408
<b>Expiration:</b>	13-May-2021

Component	Part Number	Lot Number
Bead Reagent	101535	015409
Detector Reagent	101533	015203
SBG Reagent	101534	015201
Controls	101543-101544	015501
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	1.62x10 <sup>7</sup>
Bead % Monomeric	92%
Isingle	119

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel*

11 Jun 2020

Name

Title

Signature/Date