

<b>Product Number:</b>	101537
<b>Lot Number:</b>	502723
<b>Expiration:</b>	1-Dec-2021

Component	Part Number	Lot Number
Bead Reagent	101535	103407
Detector Reagent	101533	103402
SBG Reagent	101534	103406
Control 1	101543	103502
Control 2	101544	103502
RGP Reagent	103159	N/A <sup>1</sup>

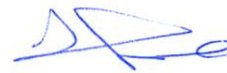
<sup>1</sup> RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.60 x10 <sup>7</sup>
Bead % Monomeric	93%
Isingle	136

Review/Approval

Junpil Park

QC Supervisor



12 Feb 2021

Name

Title

Signature/Date