

<b>Product Number:</b>	103591
<b>Lot Number:</b>	503149
<b>Expiration:</b>	16-Sep-2022

Component	Part Number	Lot Number
Bead Reagent	103581	126804
Detector Reagent	103583	126805
SBG Reagent	103585	126806
Control 1	103589	127002
Control 2	103590	127002
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	2.00 x 10 <sup>7</sup>
Bead % Monomeric	94%
Isingle	167

Review/Approval

Junpil Park

QC Supervisor

 07062024

Name

Title

Signature/Date