

<b>Product Number:</b>	103591
<b>Lot Number:</b>	502784
<b>Expiration:</b>	11-Feb-2022

Component	Part Number	Lot Number
Bead Reagent	103581	105904
Detector Reagent	103583	106016
SBG Reagent	103585	105902
Control 1	103589	106011
Control 2	103590	106011
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	1.98x10 <sup>7</sup>
Bead % Monomeric	93%
Isingle	150

Review/Approval

Junpil Park

QC Supervisor

 12 Mar 2021

Name

Title

Signature/Date