

<b>Product Number:</b>	103591
<b>Lot Number:</b>	503026
<b>Expiration:</b>	1-Jun-2022

Component	Part Number	Lot Number
Bead Reagent	103581	118104
Detector Reagent	103583	118211
SBG Reagent	103585	118212
Control 1	103589	118209
Control 2	103590	118209
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	2.02 x 10 <sup>7</sup>
Bead % Monomeric	94%
Isingle	131

Review/Approval

Junpil Park

QC Supervisor



13 Jul 2021

Name

Title

Signature/Date