

Product Number:	103591
Lot Number:	503071
Expiration:	10-Jul-2022

Component	Part Number	Lot Number
Bead Reagent	103581	119202
Detector Reagent	103583	119105
SBG Reagent	103585	119106
Control 1	103589	119404
Control 2	103590	120312
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	1.97 x 10 <sup>7</sup>
Bead % Monomeric	94%
Isingle	162

Review/Approval

Junpil Park

QC Supervisor

 30 Jul 2021

Name

Title

Signature/Date