

Product Number:	103591
Lot Number:	502771
Expiration:	11-Jan-2022

Component	Part Number	Lot Number
Bead Reagent	103581	105009
Detector Reagent	103583	105202
SBG Reagent	103585	105301
Control 1	103589	105204
Control 2	103590	105204
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	1.97x10 <sup>7</sup>
Bead % Monomeric	93%
Isingle	153

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel* 04Mar2021

Name

Title

Signature/Date