

Product Number:	103591
Lot Number:	503865
Expiration:	6-Jun-2024

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	103581	315708	315708A
Detector Reagent	103583	315805	315805A
SBG Reagent	103585	317111	317111A
Control 1	103589	316602	316602A
Control 2	103590	316602	316602A
RGP Reagent	103159	N/A <sup>1</sup>	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	92%
Isingle	145

Review/Approval

Sonia Khillou      QC      Sonia Khillou 22 JUN 2023

Name

Title

Signature/Date