

Product Number:	103591
Lot Number:	503764
Expiration:	4-Jan-2024

Component	Part Number	Lot Number
Bead Reagent	103581	301802
Detector Reagent	103583	302003
SBG Reagent	103585	302004
Control 1	103589	302304
Control 2	103590	302304
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	1.95 x 10 <sup>7</sup>
Bead % Monomeric	92%
Isingle	119

Review/Approval

*Sonia Khillan*  
Name

QC

Title

*Jan 12/16* 01FEB2023  
Signature/Date