

Product Number:	103591
Lot Number:	503324
Expiration:	5-Sep-2022

Component	Part Number	Lot Number
Bead Reagent	103581	124901
Detector Reagent	103583	125012
SBG Reagent	103585	124803
Control 1	103589	125001
Control 2	103590	125001
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	1.98 x 10 <sup>7</sup>
Bead % Monomeric	92%
Isingle	170

Review/Approval	<u>Shivani Goel</u>	QC	<u>Shivani Goel 26 Jan 2022</u>
	Name	Title	Signature/Date