

<b>Product Number:</b>	101537
<b>Lot Number:</b>	502484
<b>Expiration:</b>	23-Jun-2021

Component	Part Number	Lot Number
<b>Bead Reagent</b>	101535	021107
<b>Detector Reagent</b>	101533	021004
<b>SBG Reagent</b>	101534	021005
<b>Control 1</b>	101543	021601
<b>Control 2</b>	101544	021601
<b>RGP Reagent</b>	103159	N/A <sup>1</sup>

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

<b>Bead Count (beads/mL)</b>	1.58x10 <sup>7</sup>
<b>Bead % Monomeric</b>	90%
<b>Isingle</b>	144

<b>Review/Approval</b>	Shivani Goel	QC Manager	Shivani Goel	11 Aug 2020
	<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>	