

<b>Product Number:</b>	101537
<b>Lot Number:</b>	501847
<b>Expiration:</b>	17-Dec-2019

Component	Part Number	Lot Number
Bead Reagent	101535	835201
Detector Reagent	101533	835105
SBG Reagent	101534	835104
Controls	101543-101544	911601
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	1.61 x 10 <sup>7</sup>
Bead % Monomeric	93%
Isingle	109

Review/Approval	Patrizia Stadler	Assoc. Director QC	<i>P. Stadler</i>	07 Jun 2019
	<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>	