

Product Number:	101537
Lot Number:	501855
Expiration:	18-Apr-2020

Component	Part Number	Lot Number
Bead Reagent	101535	915801
Detector Reagent	101533	910806
SBG Reagent	101534	910807
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)	1.60 x 10 <sup>7</sup>
Bead % Monomeric	93%
Isingle	132

Review/Approval	Patrizia Stadler	Assoc. Director QC	<i>P. Stadler</i>	<i>14 Jun 2019</i>
	<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>	