

Product Number:	101537
Lot Number:	501995
Expiration:	19-Aug-2020

Component	Part Number	Lot Number
Bead Reagent	101535	923205
Detector Reagent	101533	923101
SBG Reagent	101534	923102
Controls	101543-101544	923301
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	1.62x10 <sup>7</sup>
Bead % Monomeric	93%
Isingle	142

Review/Approval

Patrizia Stadler	Assoc. Director QC	<i>p. Stadler</i>	<i>06 Sep 2019</i>
Name	Title	Signature/Date	