

Product Number:	101537
Lot Number:	501383
Expiration:	26-Jan-2019

Component	Item Number	Lot Number
Bead Reagent	101535	823305
Detector Reagent	101533	823207
SBG Reagent	101534	822805
Controls	101543-101544	821801 820701
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	1.59 x 10 <sup>7</sup>
Bead % Monomeric	93%

Review/Approval

Sivasankari Murugar QC Supervisor  
Name Title

  
Signature/Date 27 Aug 2018