

<b>Product Number:</b>	101537
<b>Lot Number:</b>	501431
<b>Expiration:</b>	16-Feb-2019

Component	Item Number	Lot Number
Bead Reagent	101535	826701
Detector Reagent	101533	826201
SBG Reagent	101534	822809
Controls	101543-101544	827101
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	1.62 x 10 <sup>7</sup>
Bead % Monomeric	93%

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

  
Signature/Date