

<b>Product Number:</b>	101537
<b>Lot Number:</b>	501563
<b>Expiration:</b>	20-Jun-2019

Component	Item Number	Lot Number
Bead Reagent	101535	835201
Detector Reagent	101533	835105
SBG Reagent	101534	835104
Controls	101543-101544	835401
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%

Review/Approval

Sivasankari Murugan

QC Supervisor

*Sivasankari Murugan* 28DEC2018

Name

Title

Signature/Date