

<b>Product Number:</b>	103187
<b>Lot Number:</b>	501479
<b>Expiration:</b>	25-Apr-2019

Component	Item Number	Lot Number
Rainbow Beads	101069	MF-030518-04
Low AEB Beads	103188	823202
High AEB Beads	101070	823203
RGP Reagent	103159	N/A <sup>1</sup>

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

**Rainbow Beads**

<b>Bead Count (beads/mL)</b>	2.39 x 10 <sup>6</sup>
<b>Bead % Monomeric</b>	91%

**Low AEB Beads**

<b>Bead Count (beads/mL)</b>	2.40 x 10 <sup>6</sup>
<b>Bead % Monomeric</b>	90%


**High AEB Beads**

<b>Bead Count (beads/mL)</b>	2.39 x 10 <sup>6</sup>
<b>Bead % Monomeric</b>	92%

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

 31 OCT 2018  
Signature/Date