

Product Number:	103187
Lot Number:	501487
Expiration:	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**

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

**Low AEB Beads**

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

**High AEB Beads**

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

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

 12/NOV/2018  
Signature/Date