

<b>Product Number:</b>	103187
<b>Lot Number:</b>	501839
<b>Expiration:</b>	7-Sep-19

Component	Item Number	Lot Number
Rainbow Beads	101069	902403
Low AEB Beads	103188	834604
High AEB Beads	101070	835503
RGP Reagent	103159	N/A <sup>1</sup>

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

**Rainbow Beads**

Bead Count (beads/mL)	2.44 x 10 <sup>6</sup>
Bead % Monomeric	86%

**Low AEB Beads**

Bead Count (beads/mL)	2.52 x 10 <sup>6</sup>
Bead % Monomeric	86%

**High AEB Beads**

Bead Count (beads/mL)	2.37 x 10 <sup>6</sup>
Bead % Monomeric	87%

**Review/Approval**

Patrizia Stadler Assoc. Director, QC

*P. Stadler* 21 May 2019

Name

Title

Signature/Date