

Product Number: 103187
Lot Number: 501770
Expiration: 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 ¹

¹ RGP is not Kit Lot Specific

Rainbow Beads

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

Low AEB Beads

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

High AEB Beads

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

Review/Approval

Patrizia Stadler Assoc. Director, QC

P. Stadler

16 Apr 2019

Name

Title

Signature/Date