

Product Number:	103187
Lot Number:	501870
Expiration:	12-Dec-19

Component	Item Number	Lot Number
Rainbow Beads	101069	902403
Low AEB Beads	103188	834604
High AEB Beads	101070	914202
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Rainbow Beads

Bead Count (beads/mL)	2.48 x 10 ⁶
Bead % Monomeric	92%

Low AEB Beads

Bead Count (beads/mL)	2.48 x 10 ⁶
Bead % Monomeric	85%

High AEB Beads

Bead Count (beads/mL)	2.38 x 10 ⁶
Bead % Monomeric	92%

Review/Approval

Patrizia Stadler Assoc. Director, QC

Name

Title

Signature/Date

P. Stadler

20 JUN 2019