

Product Number:	103187
Lot Number:	501889
Expiration:	01-Jan-2020

Component	Item Number	Lot Number
Rainbow Beads	101069	917616
Low AEB Beads	103188	916308
High AEB Beads	101070	917015
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	93%

Low AEB Beads

Bead Count (beads/mL)	2.51 x 10 ⁶
Bead % Monomeric	92%

High AEB Beads

Bead Count (beads/mL)	2.34 x 10 ⁶
Bead % Monomeric	92%

Review/Approval

Patrizia Stadler Assoc. Director, QC

P. Stadler

09 JUL 2019

Name

Title

Signature/Date