

Product Number:	103187
Lot Number:	501579
Expiration:	8-Jul-19

Component	Item Number	Lot Number
Rainbow Beads	101069	MF-060718-04
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)	2.45 x 10 ⁶
Bead % Monomeric	89.0%

Low AEB Beads

Bead Count (beads/mL)	2.40 x 10 ⁶
Bead % Monomeric	89.0%

High AEB Beads

Bead Count (beads/mL)	2.51 x 10 ⁶
Bead % Monomeric	88.9%

Review/Approval

Patrizia Stadler

QC Associate
Director

p. stadler

1/11/2019

Name

Title

Signature/Date