

Product Number:	103187
Lot Number:	501361
Expiration:	6-Feb-2019

Component	Item Number	Lot Number
Rainbow Beads	101069	MF-030518-04
Low AEB Beads	103188	MF-061318-05
High AEB Beads	101070	MF-061318-01
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Rainbow Beads

Bead Count (beads/mL)	2.35 x 10 ⁶
Bead % Monomeric	92%

Low AEB Beads

Bead Count (beads/mL)	2.48 x 10 ⁶
Bead % Monomeric	93%

High AEB Beads

Bead Count (beads/mL)	2.42 x 10 ⁶
Bead % Monomeric	92%

James Drumme
Senior Director, Manufacturing Operations
On behalf of:

Sivasankari Murugan QC Supervisor
Name Title

 Aug 10, 2018
Signature/Date

Review/Approval