

Product Number:	101537
Lot Number:	501106
Expiration:	26-Jul-2018

Component	Item Number	Lot Number
Bead Reagent	101535	80290
Detector Reagent	101533	80291
SBG Reagent	101534	80266
Controls	101543-101544	80310
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.61 x 10 ⁷
Bead % Monomeric	92.8%

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 07 FEB 2018
Signature/Date