

Product Number:	101537
Lot Number:	501346
Expiration:	20-Jan-2019

Component	Item Number	Lot Number
Bead Reagent	101535	821405
Detector Reagent	101533	820102
SBG Reagent	101534	820402
Controls	101543-101544	821801/820701
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.60 x 10 ⁷
Bead % Monomeric	92.9%

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 08 Aug 2018
Signature/Date