

Product Number:	101537
Lot Number:	501257
Expiration:	25-Oct-2018

Component	Item Number	Lot Number
Bead Reagent	101535	81370
Detector Reagent	101533	81154
SBG Reagent	101534	81155
Controls	101543-101544	81380
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.58 x 10 ⁷
Bead % Monomeric	93.0%

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 08 Jun 2018
Signature/Date