

Product Number:	101537
Lot Number:	501480
Expiration:	19-Apr-2018

Component	Item Number	Lot Number
Bead Reagent	101535	831210
Detector Reagent	101533	829804
SBG Reagent	101534	829210
Controls	101543-101544	831301
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.59 x 10 ⁷
Bead % Monomeric	93%

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 15 Nov 2018
Signature/Date