

Product Number:	101537
Lot Number:	501171
Expiration:	27-Aug-2018

Component	Item Number	Lot Number
Bead Reagent	101535	80650
Detector Reagent	101533	80590
SBG Reagent	101534	80580
Controls	101543-101544	80720
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.61 x 10 ⁷
Bead % Monomeric	93.0%

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 16 APR 2018
Signature/Date