


Product Number:	101537
Lot Number:	502003
Expiration:	18-Jul-2020

Component	Part Number	Lot Number
Bead Reagent	101535	919907
Detector Reagent	101533	919902
SBG Reagent	101534	919903
Controls	101543-101544	920003
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.59x10 ⁷
Bead % Monomeric	90%
Isingle	130

Review/Approval

Daniel Ritchie	Senior QC Analyst	
Name	Title	Signature/Date

*Signing on behalf of Patrizia Stadler