

Product Number:	101537
Lot Number:	502002
Expiration:	18-Jul-2020

Component	Part Number	Lot Number
Bead Reagent	101535	921306
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)	1.60x10 ⁷
Bead % Monomeric	92%
Isingle	124

Review/Approval

Daniel Ritchie

Senior QC Analyst

Patrizia Stadler / 26 Aug 2019

Name

Title

Signature/Date

*Signing on behalf of Patrizia Stadler