

Product Number:	101537
Lot Number:	502604
Expiration:	2-Nov-2021

Component	Part Number	Lot Number
Bead Reagent	101535	030714
Detector Reagent	101533	030707
SBG Reagent	101534	030704
Control 1	101543	030801
Control 2	101544	030801
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.61x10 <sup>7</sup>
Bead % Monomeric	92%
I <sub>single</sub>	130

Review/Approval

Junpil Park

QC Supervisor



11 Jan 2021

Name

Title

Signature/Date