

Product Number:	103591
Lot Number:	502840
Expiration:	30-Mar-2022

Component	Part Number	Lot Number
Bead Reagent	103581	108901
Detector Reagent	103583	108902
SBG Reagent	103585	108903
Control 1	103589	109006
Control 2	103590	109006
RGP Reagent	103159	N/A <sup>1</sup>

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

Bead Count (beads/mL)	2.04 x 10 <sup>7</sup>
Bead % Monomeric	92%
Isingle	142

Review/Approval

Junpil Park

QC Supervisor



08Apr 2021

Name

Title

Signature/Date