

Product Number:	103591
Lot Number:	502702
Expiration:	15-Dec-2021

Component	Part Number	Lot Number
Bead Reagent	103581	102005
Detector Reagent	103583	101903
SBG Reagent	103585	102205
Control 1	103589	102402
Control 2	103590	102402
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.98 x10 ⁷
Bead % Monomeric	95%
Isingle	125

Review/Approval

Junpil Park

QC Supervisor

 02 Feb 2021

Name

Title

Signature/Date