

Product Number:	103591
Lot Number:	502906
Expiration:	7-Apr-2022

Component	Part Number	Lot Number
Bead Reagent	103581	111503
Detector Reagent	103583	111504
SBG Reagent	103585	111505
Control 1	103589	111701
Control 2	103590	111701
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.98 x 10 ⁷
Bead % Monomeric	94%
Isingle	135

Review/Approval

Junpil Park

QC Supervisor

 05 May 2021

Name

Title

Signature/Date