

Product Number:	103591
Lot Number:	503128
Expiration:	12-Aug-2022

Component	Part Number	Lot Number
Bead Reagent	103581	124901
Detector Reagent	103583	125012
SBG Reagent	103585	124803
Control 1	103589	125001
Control 2	103590	125001
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.98 x 10 ⁷
Bead % Monomeric	92%
Isingle	170

Review/Approval

Junpil Park

QC Supervisor



15 Sep 2021

Name

Title

Signature/Date