

Product Number:	103591
Lot Number:	502989
Expiration:	10-May-2022

Component	Part Number	Lot Number
Bead Reagent	103581	116304
Detector Reagent	103583	116305
SBG Reagent	103585	116306
Control 1	103589	116910
Control 2	103590	116910
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	2.02 x 10 ⁷
Bead % Monomeric	93%
Isingle	149

Review/Approval

Junpil Park

QC Supervisor

 25 Jun 2021

Name

Title

Signature/Date