

Product Number:	103591
Lot Number:	503936
Expiration:	13-Sep-2024

Component	Part Number	Lot Number
Bead Reagent	103581	325604
Detector Reagent	103583	325705
SBG Reagent	103585	325806
Control 1	103589	325808
Control 2	103590	325808
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	2.02 x 10 ⁷
Bead % Monomeric	92%
Isingle	119

Review/Approval Sonia Khillan QC [Signature] 22SEP2023
 Name Title Signature/Date