

Product Number:	103591
Lot Number:	504051
Expiration:	9-Nov-2024

Component	Part Number	Lot Number
Bead Reagent	103581	331305
Detector Reagent	103583	331705
SBG Reagent	103585	331412
Control 1	103589	331408
Control 2	103590	331408
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.96 x 10 ⁷
Bead % Monomeric	93%
Isingle	124

Review/Approval Sonia Khillan QC Sonia Khillan 01DEC2023

Name Title Signature/Date