

Product Number:	103591
Lot Number:	503723
Expiration:	27-Oct-2023

Component	Part Number	Lot Number
Bead Reagent	103581	230412
Detector Reagent	103583	230502
SBG Reagent	103585	230511
Control 1	103589	230503
Control 2	103590	230503
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	155

Review/Approval Matt Koubouris QC Supervisor [Signature] 11 Nov 2022

Name

Title

Signature/Date