

Product Number:	103591
Lot Number:	503672
Expiration:	3-Aug-2023

Component	Part Number	Lot Number
Bead Reagent	103581	222306
Detector Reagent	103583	222402
SBG Reagent	103585	222403
Control 1	103589	226204
Control 2	103590	226204
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	2.00 x 10 ⁷
Bead % Monomeric	91%
Isingle	146

Review/Approval

Sonia Khillan

QC

Sonia Khillan / 27SEP2022

Name

Title

Signature/Date