

Product Number:	103591
Lot Number:	504002
Expiration:	13-Oct-2024

Component	Part Number	Lot Number
Bead Reagent	103581	328605
Detector Reagent	103583	328907
SBG Reagent	103585	328908
Control 1	103589	329001
Control 2	103590	329001
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.95 x 10 ⁷
Bead % Monomeric	93%
Isingle	135

Review/Approval

Sonia Khillan

QC

Sonia Khillan 25 OCT 2023

Name

Title

Signature/Date