

Product Number:	103591
Lot Number:	503841
Expiration:	28-Mar-2024

Component	Part Number	Lot Number
Bead Reagent	103581	308709
Detector Reagent	103583	308706
SBG Reagent	103585	308707
Control 1	103589	308710
Control 2	103590	308710
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.95 x 10 ⁷
Bead % Monomeric	91%
Isingle	126

Review/Approval

Sonia Khilken

QC

Sonia Khilken

11 APR 2023

Name

Title

Signature/Date