

Product Number:	103591
Lot Number:	504111
Expiration:	15-Feb-2025

Component	Part Number	Lot Number
Bead Reagent	103581	404011
Detector Reagent	103583	404313
SBG Reagent	103585	404604
Control 1	103589	409202
Control 2	103590	409202
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Bead Count (beads/mL)	1.97 x 10 ⁷
Bead % Monomeric	93%
Isingle	146

Review/Approval

Cody Carlson

QC Manager

Cody Carlson 05 APR 2024

Name

Title

Signature/Date