

Product Number:	103591
Lot Number:	504234
Expiration:	10-Oct-2025

Component	Part Number	Lot Number
Bead Reagent	103581	410903
Detector Reagent	103583	411001
SBG Reagent	103585	411002
Control 1	103589	411501
Control 2	103590	411501
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Bead Count (beads/mL)	2.04 x 10 ⁷
Bead % Monomeric	92%
Isingle	118


Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 09MAY2024

Signature/Date