

Product Number:	103591
Lot Number:	503070
Expiration:	24-Jul-2022

Component	Part Number	Lot Number
Bead Reagent	103581	120506
Detector Reagent	103583	120601
SBG Reagent	103585	120602
Control 1	103589	122005
Control 2	103590	122005
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.96 x 10 ⁷
Bead % Monomeric	95%
Isingle	146

Review/Approval

Wenkui Tsui

QC



18 Aug 2021

Name

Title

Signature/Date