


Product Number:	103591
Lot Number:	502946
Expiration:	30-Apr-2022

Component	Part Number	Lot Number
Bead Reagent	103581	114201
Detector Reagent	103583	114302
SBG Reagent	103585	114303
Control 1	103589	114404
Control 2	103590	114404
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	2.03 x 10 ⁷
Bead % Monomeric	94%
Isingle	153

Review/Approval	Junpil Park	QC Supervisor	 02 Jun 2021
	Name	Title	Signature/Date