


Product Number:	103591
Lot Number:	503476
Expiration:	16-Feb-2023

Component	Part Number	Lot Number
Bead Reagent	103581	213605
Detector Reagent	103583	213601
SBG Reagent	103585	213602
Control 1	103589	213701
Control 2	103590	213701
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.96 x 10 ⁷
Bead % Monomeric	92%
Isingle	151

Review/Approval	Matthew Smart	QC Director	 03 Jun 2022
	Name	Title	Signature/Date