

Product Number:	103591
Lot Number:	503201
Expiration:	30-Oct-2022

Component	Part Number	Lot Number
Bead Reagent	103581	130306
Detector Reagent	103583	130303
SBG Reagent	103585	130307
Control 1	103589	134114
Control 2	103590	134114
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	2.00 x 10 ⁷
Bead % Monomeric	94%
I _{single}	131

Review/Approval	<u>Shivani Goel</u>	QC	<u>Shivani Goel 30 Dec 2021</u>
	Name	Title	Signature/Date