

Product Number:	101537
Lot Number:	502604
Expiration:	22-Oct-2021

Component	Part Number	Lot Number
Bead Reagent	101535	030714
Detector Reagent	101533	030707
SBG Reagent	101534	030704
Control 1	101543	030801
Control 2	101544	030801
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.61x10 ⁷
Bead % Monomeric	92%
Isingle	130

Review/Approval Shivani Goel QC Manager
Name Title

Shivani Goel 12NOV2020
Signature/Date