

Component	Part Number	Lot Number
Bead Reagent	101535	919907
Detector Reagent	101533	919902
SBG Reagent	101534	919903
Controls	101543-101544	920003
RGP Reagent	103159	N/A ¹

Bead Count (beads/mL)	1.59×10^7
Bead % Monomeric	90%
I _{single}	130