

Product Number:	101537
Lot Number:	501878
Expiration:	18-Jun-2020

Component	Part Number	Lot Number
Bead Reagent	101535	917603
Detector Reagent	101533	917103
SBG Reagent	101534	917006
Controls	101543-101544	917701
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.59 x 10 ⁷
Bead % Monomeric	92%
Isingle	125

Review/Approval	Patrizia Stadler	Assoc. Director	<i>P. Stadler</i>	11 JUL 2019
		QC		
	Name	Title	Signature/Date	