

Product Number:	101537
Lot Number:	502050
Expiration:	20-Sep-2020

Component	Part Number	Lot Number
Bead Reagent	101535	926802
Detector Reagent	101533	926403
SBG Reagent	101534	926302
Controls	101543-101544	926901
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.58x10 ⁷
Bead % Monomeric	93%
Isingle	155

Review/Approval	Patrizia Stadler	Assoc. Director QC	<i>P. Stodler</i>	08 OCT 2019
	Name	Title	Signature/Date	