

Product Number:	101537
Lot Number:	501751
Expiration:	14-Mar-2020

Component	Item Number	Lot Number
Bead Reagent	101535	909114
Detector Reagent	101533	907904
SBG Reagent	101534	907304
Controls	101543-101544	910201
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.61 x 10 ⁷
Bead % Monomeric	93%
Isingle	154

Review/Approval	Patrizia stadler	QC Assoc. Director	<i>P Stadler</i>	17 Apr 2019
	Name	Title	Signature/Date	