

Product Number:	101537
Lot Number:	502302
Expiration:	13-Mar-2021

Component	Part Number	Lot Number
Bead Reagent	101535	009101
Detector Reagent	101533	008505
SBG Reagent	101534	008302
Controls	101543-101544	009202
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.61x10 ⁷
Bead % Monomeric	93%
Isingle	136

Review/Approval

	Assoc. Director		
Patrizia Stadler	QC	<i>P. Stadler</i>	08 Apr 2020
Name	Title	Signature/Date	