

Product Number:	101537
Lot Number:	502124
Expiration:	25-Aug-2020

Component	Part Number	Lot Number
Bead Reagent	101535	932301
Detector Reagent	101533	932202
SBG Reagent	101534	932201
Controls	101543-101544	932401
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.60x10 ⁷
Bead % Monomeric	93%
Isingle	178

Review/Approval

Patrizia Stadler	Assoc. Director QC	<i>P. Stadler</i>	27 Nov 2019
Name	Title	Signature/Date	