

Product Number:	101537
Lot Number:	501732
Expiration:	14-Mar-2020

Component	Item Number	Lot Number
Bead Reagent	101535	908005
Detector Reagent	101533	907904
SBG Reagent	101534	907304
Controls	101543-101544	908101
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.59x10 ⁷
Bead % Monomeric	93%

Review/Approval

Patrizia stadler QC Assoc. Director

P. Stadler 28 Nov 2019

Name

Title

Signature/Date