

Product Number:	101537
Lot Number:	501668
Expiration:	8-Feb-2020

Component	Item Number	Lot Number
Bead Reagent	101535	905006
Detector Reagent	101533	904503
SBG Reagent	101534	903904
Controls	101543-101544	905101
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.62 x 10 ⁷
Bead % Monomeric	99%

Review/Approval

Patrizia stadler

QC Assoc. Director

P. Stadler 26 Feb 2019

Name

Title

Signature/Date