

Product Number:	103591
Lot Number:	502316
Expiration:	5-Apr-2021

Component	Part Number	Lot Number
Bead Reagent	103581	023108
Detector Reagent	103583	009601
SBG Reagent	103585	009602
Control 1	103589	025908
Control 2	103590	025908
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	2.00x10 ⁷
Bead % Monomeric	91%
Isingle	168

Review/Approval Patrizia Stadler Assoc. Director, QC *P. Stadler* 29 Sep 2020
 Name Title Signature/Date