

Product Number:	101537
Lot Number:	502204
Expiration:	27-Dec-2020

Component	Part Number	Lot Number
Bead Reagent	101535	002113
Detector Reagent	101533	002201
SBG Reagent	101534	002102
Controls	101543-101544	002301
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Bead Count (beads/mL)	1.58x10 ⁷
Bead % Monomeric	93%
Isingle	135

Review/Approval	Assoc. Director		
	QC	<i>P. Stadler</i>	<i>3 Feb 2020</i>
	<u>Patrizia Stadler</u>	<u></u>	<u></u>
	Name	Title	Signature/Date