

Product Number:	101354
Lot Number:	502307
Expiration:	18-Dec-2020

Component	Item Number	Lot Number
Beads	103207	003610
Detector/Sample Diluent	101359	007402
Biotinylation Reaction Buffer	101358	935220
Bead Conjugation Buffer	101357	001506
Bead Blocking Buffer	101356	935219
Bead Wash Buffer	101355	008007
Bead Diluent Buffer	100458	006215

Bead Count (beads/mL)	2.14x10 ⁹
--------------------------	----------------------

Review/Approval

Patrizia Stadler	QC Assoc. Director	<i>p. stadler</i> 02Apr 2020
Name	Title	Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.