

Product Number:	101354
Lot Number:	502182
Expiration:	1-Aug-2020

Component	Item Number	Lot Number
Beads	103207	912114
Detector/Sample Diluent	101359	932502
Biotinylation Reaction Buffer	101358	935207
Bead Conjugation Buffer	101357	935108
Bead Blocking Buffer	101356	935209
Bead Wash Buffer	101355	932505
Bead Diluent Buffer	100458	921319

Bead Count (beads/mL)	3.27x10 ⁹
--------------------------	----------------------

Review/Approval

Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i> 31 Dec 2019
Name	Title	Signature/Date

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