

Product Number:	101354
Lot Number:	502367
Expiration:	9-Oct-2020

Component	Item Number	Lot Number
Beads	103612	005516
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	928215

Bead Count (beads/mL)	3.07x10 ⁹
--------------------------	----------------------

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 11 May 2020

Name

Title

Signature/Date

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