

Product Number:	101354
Lot Number:	502471
Expiration:	18-Dec-2020

Component	Item Number	Lot Number
Beads	103612	005518
Detector/Sample Diluent	101359	015806
Biotinylation Reaction Buffer	101358	015108
Bead Conjugation Buffer	101357	011706
Bead Blocking Buffer	101356	935219
Bead Wash Buffer	101355	015106
Bead Diluent Buffer	100458	007710

Bead Count (beads/mL)	3.30x10 ⁹
--------------------------	----------------------

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 20/01/2020
Signature/Date

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