

Product Number:	101354
Lot Number:	502494
Expiration:	24-May-2021

Component	Item Number	Lot Number
Beads	103612	005518
Detector/Sample Diluent	101359	019711
Biotinylation Reaction Buffer	101358	015108
Bead Conjugation Buffer	101357	015107
Bead Blocking Buffer	101356	019812
Bead Wash Buffer	101355	015106
Bead Diluent Buffer	100458	014512

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

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 19 Aug 2020

Name

Title

Signature/Date

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