

Product Number:	101354
Lot Number:	502443
Expiration:	18-Dec-2020

Component	Item Number	Lot Number
Beads	103612	015310
Detector/Sample Diluent	101359	015806
Biotinylation Reaction Buffer	101358	015108
Bead Conjugation Buffer	101357	011706
Bead Blocking Buffer	101356	935219
Bead Wash Buffer	101355	008007
Bead Diluent Buffer	100458	007710

Bead Count (beads/mL)	3.36x10 ⁹
--------------------------	----------------------

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 29 Jun 2020

Name

Title

Signature/Date

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