

Product Number:	101354
Lot Number:	502526
Expiration:	30-May-2021

Component	Item Number	Lot Number
Beads	103612	015313
Detector/Sample Diluent	101359	019711
Biotinylation Reaction Buffer	101358	023711
Bead Conjugation Buffer	101357	015107
Bead Blocking Buffer	101356	019812
Bead Wash Buffer	101355	015106
Bead Diluent Buffer	100458	022412

Bead Count (beads/mL)	3.24x10 ⁹
--------------------------	----------------------

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 10 Sep 2020
Signature/Date

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